

Introduction to

# Cardiovascular Disease



Of course, we've taken liberties on the cover: jellyfish don't have hearts. The surface of the bell-shaped body is used for gas exchange, so even respiratory and vascular systems are unneeded. Digestion involves only a simple cavity, lined with a gastrodermis, for both food intake and waste expulsion. Some jellyfish move by contracting muscles that line the bell, expelling water and propelling the body forward. For these jellies, such movement is largely for repositioning in the water column, guided by rudimentary sensors for light and gravity. However, the box jellyfish, of the class *Cubozoa*, is another beast altogether. These hunters have an umbrella-shaped body with true eyes, complete with retinas, corneas, and lenses. Improved vision, combined with neural and structural modifications of the body, allows rapid, directional (and, presumably, intentional) movement. Perhaps most importantly, box jellyfish tentacles are studded with cells containing barbed, pressurized nematocysts that inject venom into the victim. The venom of the most dangerous box jelly, *Chironex fleckeri*, has dermonecrotic, myotoxic, cardiotoxic, hemodynamic, and hemolytic effects. Envenomation, particularly in smaller children, can be lethal. More commonly, symptoms include severe localized pain, marked redness of the skin, and collapse due to respiratory failure or cardiac arrest.

Compared to an encounter with a *Chironex*, cardiovascular disease might seem relatively mild. However, statistics from the American Heart Association indicate that in 2006, 6,161,000 Americans were discharged from short hospital stays with a first listed diagnosis of cardiovascular disease. Roughly 40% of men and women in the 40 to 59 year old age group have cardiovascular disease. This increases to 73% percent of men and women in the 60-79 year old age group. For both men and women, black, white, or Asian, cardiovascular diseases caused more deaths than cancer in 2006. More than 800,000 people died from cardiovascular diseases in the United States in 2006, while some 70 people have died from *Chironex* poisoning in recorded history.

The statistics on deaths due to cardiovascular disease are startling, but the overall cost, in terms of treatment and diminished quality of life, must be staggering. Cayman is dedicated to helping make research possible, by making quality reagents available at affordable prices. Our product line extends well beyond the contents of this topical catalog. Please look at our website, caymanchem.com, or call us to find out how we can work together to understand cardiovascular disease.



Cardiovascular Disease, Pulmonary Hypertension, and Prostacyclins

Epigenetics: The New Frontier for Vascular Diseases

You Say 'Thrombosis', I Say 'Thromboxane'

table of contents

Think of Cayman as Epigenetics Central

eNOS, Vascular Endothelial Health, and CVD



Orders are accepted by telephone, fax, mail, e-mail, or via the Cayman Chemical website. We will accept telephone orders Monday through Friday from 8 AM to 6 PM EST. All orders received by 1 PM EST will be shipped the same day if stock is available (Monday through Thursday only). Confirming purchase orders must be clearly marked as such to avoid possibility of duplication.

ordering information

In most instances we ship FedEx Standard Overnight Delivery (not available to all locations), with delivery by 3:30 PM of the next business day. Product availability may vary. Local delivery is available for the Ann Arbor area only. Other shipping options will be considered upon request, but can be granted only under conditions that will ensure the quality of the product. Freight is prepaid and added to the invoice. Please inquire at the time of order for an estimate of the freight charges. If you wish us to ship collect, please supply a valid account number when ordering. Please address all orders to:

Toll-free Phone: (800) 364-9897 Cayman Chemical Company Fax: (734) 971-3640 1180 E. Ellsworth Road Ann Arbor, MI 48108 USA E-mail: custserv@caymanchem.com Phone: (734) 971-3335 www.caymanchem.com

#### Include the following information with your order:

- 1. Catalog number, description, size, and quantity desired.
- 2. Complete shipping address. (Delivery is not available to post office box numbers.)
- 3. A complete billing address
- 4. A purchase order number or major credit card (Visa, MasterCard, or American Express), account number, and expiration date. 5. Name of the end user.

- 1. U.S. funds only, drawn on a U.S. bank 2. Net 30 days.
- 3. F.O.B. Ann Arbor, Michigan, U.S.A.
- 4. Bank fees and wire transfer fees are not to be deducted from the invoice amount.

All prices listed are in U.S. dollars. The prices in this catalog are effective as of May 1, 2011. Prices are subject to change without notice.

Products cannot be returned without prior authorization from Cayman Chemical Company. Please contact our Customer Service Department for return shipping instructions. Custom orders and radioactive material cannot be accepted for return credit if due to a

#### Technical Assistance

Technical assistance is available from 8 AM to 5:30 PM EST. If inquiring about a purchased product, please provide the catalog number, lot number, and date of purchase to our technical staff so they may answer your questions quickly. Technical assistance may be reached toll free at 888-526-5351, via e-mail at techserv@caymanchem.com, or on the web at www.caymanchem.com/techserv.

#### **Use of Research Products**

The products in this catalog are not for human or veterinary disease diagnosis or therapeutic drug use. They should be used only by technically qualified individuals or those under their direct supervision. Any individual working directly with these products should have free access to the applicable Material Safety Data Sheet (MSDS) and should read and understand it completely prior to use. Please contact our Customer Service Department or visit the specific product page on our website if you require additional copies of any MSDS.

The end-user assumes full responsibility for appropriate licensing and/or non-infringement for any proprietary claim or patent.

NOTE: For Laboratory Research Use Only. Not for human or veterinary diagnostic or therapeutic use.

DTT DGAT DIO DNA DMSO EDTA EET EGFR EIA ELISA EP EPA ER FXR GC GCS GPCR GSK GST HDL HIF HMG HMT HRP HSP ICC IF IHC IL bbreviations

NOS oxLDL PAF-AH PBS PE PG PGES PGIS PI3K PKA PL PP PRMT PRP PRMT PRP RNA RTK SAME SEH SGC SIP SIRT STEBP

CK CNS CoA COX CREB CYP DTNB ,5'-Dithio-bis-(2-nitrobenzoic acid); Platelet-activating Factor Platelet-activatingFactor Acetylhydrolase

helial Growth Factor

caymanchem.com

Cayman Chemical

# Cardiovascular Disease, Pulmonary Hypertension, and Prostacyclins

The term 'cardiovascular disease' can be confusing. Commonly, it centers on atherosclerosis, focusing on cholesterol levels, plaque formation, and the narrowing and hardening of the arteries. From here, it is necessarily linked to the other diseases that lead to and result from these changes, including hypercholesterolemia, diabetes, hypertension, and stroke. This broadens the scope of the relevant pathophysiology tremendously. Atherosclerosis can be further linked with metabolic syndrome, which is not so much a disease as a group of risk factors (obesity, insulin resistance, aging, genetic predisposition) that increase the risk for coronary artery disease, stroke, and type 2 diabetes. Cardiovascular disease, in this broad sense then, is not so much a single disease as an extensive web of interconnected diseases.

A different approach is given by the International Classification of Diseases (ICD), maintained by the World Health Organization. Originally called the International List of Causes of Death, it was created in the 1850s to standardize the classification of causes of morbidity worldwide. Now in its tenth revision, the ICD codes over 14,000 diseases, symptoms, complaints, social circumstances, and external causes of injury or diseases. These are divided into 22 broad categories, including the diseases of the circulatory system. Circulatory system diseases span several hundred subcategories, ranging from ischemic, rheumatic, and pulmonary heart diseases to diseases of veins and arteries to subcerebral hemorrhage and stroke. In this system of classification, several types of atherosclerosis are categorized within the Diseases of Arteries, Arterioles, and Capillaries. According to ICD-10, then, the diseases of the cardiovascular system can be neatly divided into numerous diverse and relatively distinct entities.

#### **Pulmonary Hypertension**

Hypertension is an important topic under both approaches. Hypertension is simply the technical term for high blood pressure. Blood pressure is affected, physiologically, by such factors as water and salt content, condition of kidneys, nervous system, or blood vessels, and hormone levels. Hypertension is a risk factor for severe adverse events, such as heart attack, stroke, heart failure, kidney disease, and early death. Pulmonary hypertension is specifically an increase in blood pressure in the lung vasculature and is characterized, physiologically, by dyspnea (shortness of breath), dizziness, and fainting. The right ventricle of the heart pumps blood through the lungs, where it can receive oxygen (Figure 1). When the small vessels within the lungs become narrow, the resistance to blood flow through the lungs increases, raising pulmonary blood pressure. If the narrowing is due to non-inflammatory processes, then it is called pulmonary hypertension. Because the heart works harder to move blood through the narrowed vessels, the right ventricle may become enlarged. Persistent pulmonary hypertension inevitably leads to heart failure.

A decade or more ago, pulmonary hypertension was divided into primary (essential) pulmonary hypertension, which by definition was idiopathic, and hypertension that was secondary to, commonly, other cardiac or pulmonary disorders. More recently, pulmonary hypertension has been further classified, by the World Health Organization, into five types, including arterial, venous, hypoxic, thromboembolic, and miscellaneous. Pulmonary arterial hypertension (PAH) is characterized by abnormally high blood pressure in the pulmonary artery, the blood vessel that carries blood from the heart to the lungs (Figure 1).

#### **Lipid Mediators in Pulmonary Hypertension**

In the older literature, primary pulmonary hypertension (PPH) was reported to be commonly associated with lipid mediator dysregulation. An overabundance of thromboxanes during PPH was described as indicative of persistent platelet activation, while an underproduction of PGI2 indicated

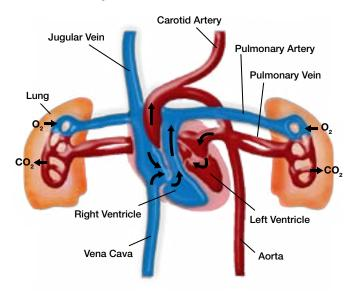


Figure 1. Deoxygenated blood returning to the heart is pumped through the right ventricle to the lungs. The smallest vessels of the lungs, which are crucial for gas exchange, are key points of resistance to blood flow and modulate pulmonary blood pressure.

an inadequate endothelial response. In addition, PPH often featured an increase in PGD<sub>2</sub>, indicating activation of tissue macrophages. In terms of the newer disease classification, these changes are featured primarily in PAH. Analogs of PGI<sub>2</sub> are effective, either alone or in combination therapy, in the treatment of pulmonary hypertension (PPH or PAH).<sup>2,3</sup> In part, this results from their ability to relax pulmonary arterial smooth muscle. Moreover, PGI<sub>2</sub> and its analogs may also suppress the overproduction of thromboxane and PGD<sub>2</sub>, since they can potently inhibit macrophage function, including their production of inflammatory mediators.<sup>4</sup>

The synthesis of PGI<sub>2</sub> from arachidonic acid is initiated by the cyclooxygenases, COX-1 and COX-2 (Figure 2). Both enzymes are membrane-bound homodimers, found predominantly on the perinuclear membranes. COX-1 is commonly described as ubiquitous and constitutively expressed. COX-2, on the other hand, is normally absent from most cell types, its expression is induced by inflammatory mediators like TNF-α and IL-1B, and both its mRNA and protein have short half-lives. Both COX isoforms produce the intermediate PGH<sub>2</sub>, which is then converted to the final prostanoid products by specific, terminal enzymes. For example, PGH<sub>2</sub> is isomerized to PGI<sub>2</sub> by PGIS (also known as CYP8A1), a singlepass membrane-bound protein that is abundant in endothelium, ovary, heart, skeletal muscle, lung, and prostate. While the expression of PGIS is constitutive, the production of PGI2 can be dramatically increased by the induced expression of COX-2, particularly in vascular cells and macrophages. PGI<sub>2</sub> is rapidly secreted from source cells, presumably through an ATP-binding cassette (ABC) transporter. PGI<sub>2</sub> activates neighboring cells which bear the specific I-prostanoid (IP) receptor, a GPCR that signals through Gas to activate adenylate cyclase (AC) and elevate intracellular cAMP levels (Figure 2). Through this pathway, PGI<sub>2</sub> relaxes vascular and bronchial smooth muscle, inhibits platelet aggregation, and suppresses the production of proinflammatory cytokines, angiogenic cytokines, and mediators of extracellular matrix remodeling (e.g., IL-1, IL-6, VEGF, TGF-β, thromboxane). In the circulation, PGI<sub>2</sub> is non-enzymatically hydrated to 6-keto PGF<sub>1 $\alpha$ </sub> (t<sub>1/2</sub> = 2-3 minutes), and then converted to the major metabolite, 2,3-dinor-6-keto PGF<sub>1 $\alpha$ </sub> ( $t_{1/2}$  = 30 minutes).

Epoprostenol is another name for PGI<sub>2</sub>. The sodium salt, epoprostenol sodium, is marketed for therapeutic use in the formulation called Flolan<sup>TM</sup>. Used in the treatment of PAH, it must be given intravenously because of its very short physiologic half-life. Carbaprostacyclin (originally called carbacyclin) is a first generation analog of PGI2 with a simple carbon-foroxygen substitution. While it is about 30 times less potent than PGI2 as an inhibitor of induced platelet aggregation, it is more chemically stable.<sup>5</sup> Surprisingly, carbaprostacyclin has been shown to be an effective activator of peroxisome proliferator-activated receptor  $\delta$  (PPAR $\delta$ ), which is involved in adipocyte regulation and modulates monocyte/macrophage function.<sup>6</sup> This suggests that PGI<sub>2</sub> may be an endogenous ligand for PPARδ, which might be particularly relevant when COX-2 is induced and elevated levels of PGI<sub>2</sub> are generated. Cicaprost is a PGI<sub>2</sub> analog that is orally active with prolonged availability in vivo, having a terminal half life in plasma of one hour. Importantly, cicaprost has been shown to strongly reduce lung and lymph node metastasis in rats, suggesting that it might be useful in cancer

Beraprost, like cicaprost, is an orally active PGI2 analog. Marketed as an ingredient in the formulations Dorner<sup>TM</sup> and Procyclin<sup>TM</sup>, it has a half-life in plasma of several hours, effectively improves pulmonary hemodynamics in patients with pulmonary hypertension, and inhibits platelet aggregation. In clinical trials, beraprost has been found to be safe and well-tolerated.9 It is now being tested for efficacy in peripheral vascular disease, Buerger's disease, arteriosclerosis obliterans, and reduction of reperfusion injury, as well as in PAH. Iloprost (ingredient of Ventavis<sup>TM</sup>), a second generation analog of PGI<sub>2</sub>, is used in aerosolized form. It binds with equal affinity to the human IP and EP<sub>1</sub> (PGE<sub>2</sub>) receptors, a K<sub>1</sub> value of 11 nM.<sup>10</sup> In whole animals, iloprost acts as a vasodilator, hypotensive, antidiuretic, and prolongs bleeding time. It has been demonstrated to be effective in the treatment of PAH and is well-tolerated in long-term usage. 11,12 Intravascular iloprost has mixed utility in the treatment of Raynaud's syndrome. 13

Treprostinil is one of the newer stable PGI<sub>2</sub> analogs and is currently used for the treatment of PAH, formulated for infusion, either subcutaneous or intravenous (Remodulin<sup>TM</sup>), and inhalation (Tyvaso<sup>TM</sup>) are available. In addition to its direct vasodilatory effects, it also potently inhibits macrophage function, preventing NF-KB activation, phagocytosis, and production of proinflammatory cytokines. 4,13

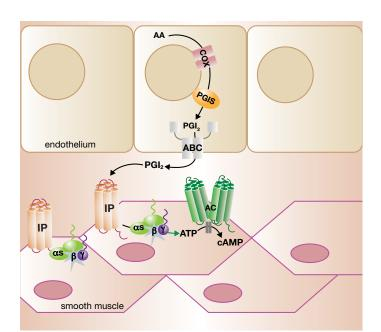


Figure 2. PGI, is synthesized from AA by the COX pathway via PGIS. Secreted PGI, activates the IP receptor, which induces cAMP synthesis from ATP by AC. In the vascular system, PGI<sub>2</sub> from the endothelium causes relaxation of vascular smooth muscle.

A common test for effectiveness of treatment is the 'six minute walk', performed in the pulmonary function facility. The test provides information on how far the patient can walk in six minutes and what happens to blood oxygen levels during exertion. With these data, treatments can be assessed and optimized. Because each of the prostacyclin analogs is distinctive in its mode of delivery, stability in plasma, and additional actions, the choice for treatment must be matched to the patient's needs.

Prostacyclin and Analogs for Research Use			
Item No.	Product Name		
16831	Cicaprost	Taken orally; half life in plasma of one hour	
18230	Beraprost	Orally active (Dorner $^{\text{TM}}$ , Procyclin $^{\text{TM}}$ ); half life of several hours in plasma	
18215	lloprost	Inhaled (Ventavis <sup>TM</sup> )	
10162	Treprostinil	Injected (Remodulin <sup>TM</sup> ) or inhaled (Tyvaso <sup>TM</sup> )	
18220	$ \begin{array}{c} \textbf{Prostaglandin I}_2 \\ \textbf{(Sodium Salt)} \end{array} \qquad \textbf{Injected (Flolan}^{\text{TM}}) $		
18210	Carbaprostacyclin	Taken orally; rapidly metabolized	
For a full product listing, please visit www.caymanchem.com			

- Christman, B.W. Chest 114(3), 205S-207S (1998)
- Galiè, N., Negro, L., and Simonneau, G. Eur. Respir. Rev. 18(113), 148-153 (2009).
- Delcroix, M., Spaas, K., and Ouarck, R. Eur. Respir, Rev. 18(117), 253-259 (2009).
- Aronoff, D.M., Peres, C.M., Serezani, C.H., et al. J. Immunol. 178, 1628-1634 (2007).
- Whittle, B.J.R. and Moncada, S. Circulation 72, 1219-1225 (1985)
- Gupta, R.A., Tan, J., Krause, W.F., et al. Proc. Natl. Acad. Sci. USA 97(24), 13275-13280 (2000).
- Hildebrand, M., Staks, T., and Nieuweboer, B. Eur. J. Clin. Pharmacol. 39(2), 149-153 (1990). Schirner, M., Kraus, C., Lichtner, R.B., et al. Prostaglandins Leukot. Essent. Fatty Acids 58(40) 311-317 (1998).
- Melian, E.B. and Goa, K.L. Drugs 62(1), 107-133 (2002).
- Ahramovitz M. Adam M. Roje Y. et al. Biochim Biophys. Acta 1483, 285-293 (2000).
- Olschewski, H., Simonneau, G., Galié, N., et al. N. Engl. J. Med. 347(5), 322-329 (2002).
- Hoeper, M.M., Schwarze, M., Ehlerding, S., et al. N. Engl. J. Med. 342, 1866-1870 (2000)
- Raychauduri, B., Malur, A., Bonfield, T.L., et al. J. Biol. Chem. 277(36), 33344-33348 (2002).

13042

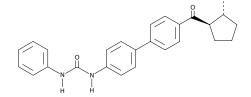
10008463

[959122-11-3] (1R,2R)-2-(4'-(3-phenylureido)biphenylcarbonyl)-Cyclopentane Carboxylic Acid)

MF:  $C_{26}H_{24}N_2O_4$  FW: 428.5 Purity:  $\geq 95\%$ A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A potent inhibitor of DGAT-1 (IC<sub>50</sub>s = 7 and 24 nM, for human and murine, respectively); confers significant weight loss within seven days and significantly reduces plasma and liver triglycerides when administered at 3 mg/kg to DIO mice

1 mg 5 mg 10 mg 25 mg

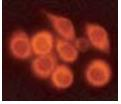


#### **ACAT-1 Polyclonal Antibody**

100028 Acyl-coenzyme A:Cholesterol Acyltransferase-1, Sterol O-Acyltransferase 1

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human ACAT-1 amino acids 6-23 • Host: rabbit • Cross Reactivity: (+) murine, rat, porcine, and human ACAT-1 • Application(s): ICC, IHC, and WB • ACAT-1 catalyzes the formation of cholesterol esters from cholesterol and long chain fatty acyl-CoA, and may play a role in the development of atherosclerosis.



• Also Available: ACAT-1 Blocking Peptide (10005090)

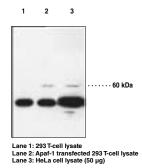
#### **ACAT-2 Polyclonal Antibody**

100027

Acyl-coenzyme A:Cholesterol Acyltransferase-2, Sterol O-Acyltransferase 2 Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human ACAT-2 amino acids 3-20 • Host: rabbit • Cross Reactivity: (+) human, murine, rat, porcine, and ovine ACAT-2 • Application(s): ICC, IHC, and WB • ACAT-2 catalyzes the formation of cholesterol esters from cholesterol and long chain fatty acyl-CoA.

500 µl



• Also Available: ACAT-2 Blocking Peptide (10005091)

#### Acenocoumarol

10010569

10007686

589451

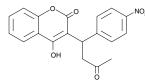
[152-72-7] Acenocoumarin, G-23350, Nicoumalone, Sinthrome, Sintron

MF:  $C_{19}H_{15}NO_6$  FW: 353.3 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An analog of the oral anticoagulant warfarin; functions as a vitamin K antagonist and has a half-life in plasma of about eight hours

50 mg 100 mg 250 mg



• Also Available: (R)-(+)-Acenocoumarol (9000336) (S)-(-)-Acenocoumarol (9000337)

#### Acetyl Podocarpic Acid Anhydride

[344327-48-6] APD

MF:  $C_{39}H_{46}O_7$  FW: 614.8 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A potent, semi-synthetic LXR agonist derived from extracts of the mayapple; induces the expression of the ABCA1 reverse cholesterol transporter to increase the efflux of cholesterol from enterocytes and thus inhibits the overall absorption of cholesterol (ED<sub>50</sub> = 1 nM)

1 mg 5 mg 10 mg 50 mg

AcSDKP EIA Kit+ N-Acetyl Ser-Asp-Lys-Pro

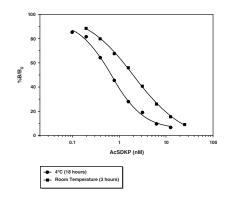
**Stability:** ≥6 months at -20°C

Limit of Detection: 1.5 pg/ml

Sensitivity: 50% B/B<sub>0</sub>: 2.0 nM after 3 hour immunological reaction • 0.5 nM after 18 hour immunological reaction • 80% B/B<sub>0</sub>: 0.2 nM after 18 hour immunological

Summary: AcSDKP is a tetrapeptide growth regulatory hormone which inhibits the proliferation of hematopoietic stem cells. The dipeptidase angiotensin converting enzyme (ACE) actively metabolizes circulating AcSDKP, giving it a brief plasma half-life of 4 to 5 minutes. ACE inhibition is a major therapeutic end point in the treatment of hypertension management. A further consequence of ACE inhibition is the accumulation of AcSDKP in plasma and urine. This accumulation may have physiological effects, which are manifested as the anemia of chronic ACE inhibitor toxicity. More commonly, plasma and urine AcSDKP levels can be used as a biomarker of ACE inhibition and an index of patient compliance with therapy. Measurement of AcSDKP in human urine or plasma can be readily accomplished by

96 wells



#### Adipose Trialyceride Lipase Polyclonal Antibody

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

catalyzing the conversion of triacylglycerols to diacylglycerols.

Summary: Antigen: human adipose triglyceride lipase (ATGL) amino acids 382-400

• Host: rabbit • Cross Reactivity: (+) human, murine, and rat ATGL • Application(s):

IHC (paraffin-embedded sections) and WB • ATGL is one of the key enzymes

involved in the mobilization of fatty acids from triglyceride stores in adipose tissue,

1 2 3 4

Lane 1:Human liver microsome (50 μg) Lane 2: HepG2 cell lysate (50 μg) Lane 3: Rat brown fat homogenate (50 μ Lane 4: Murine liver 100,000 x g pellet (50 μ

ATGL, Desnutrin, PLA, C

500 ul

10006409 [217082-58-1]

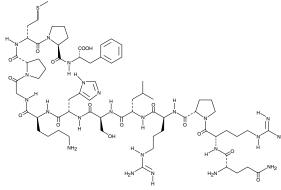
**MF:**  $C_{69}H_{111}N_{23}O_{16}S$  **FW:** 1,550.8 **Purity:**  $\geq$ 95%

Apelin-13

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: Endogenous ligand of the APJ receptor, with an EC50 value of 0.37 nM; acts primarily in the periphery and CNS, playing important roles in regulating cardiovascular function, fluid homeostasis, hypertension, and insulin sensitivity

1 mg 5 mg 10 mg 25 mg



#### • Also Available: Adipose Triglyceride Lipase Blocking Peptide (10010569)

#### AMP-Deoxynojirimycin

10010332

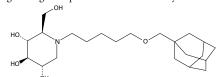
[216758-20-2] Adamantane-pentyl-dNM, AMP-dNM, N-(5-adamantane-1-ylmethoxy-pentyl)-Deoxynojirimycin

MF:  $C_{22}H_{39}NO_5$  FW: 397.6 Purity:  $\geq$ 95%

A solution in ethanol **Stability:**  $\ge 1$  year at  $-20^{\circ}$ C

Summary: A hydrophobic derivative of deoxynojirimycin that potently inhibits BGD  $(IC_{50} = 0.3 \text{ nM})$ , less potently antagonizes GCS  $(IC_{50} = 25 \text{ nM})$ , but only poorly inhibits other GCase isoforms; suppresses hapten-induced colitis, enhances insulin sensitivity, and induces SREBP-regulated gene expression and cholesterol synthesis

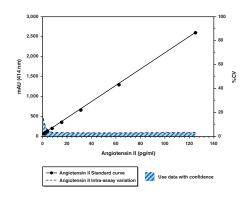


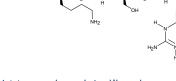


#### 589301 Angiotensin II EIA Kit

**Stability:** ≥6 months at -20°C Limit of Detection: 1.5 pg/ml

Summary: Angiotensin II is a primary reactive vasoconstrictor, the main stimulus for aldosterone release, and one of the causative factors of chronic hypertension. The active angiotensin II octapeptide is released via a tightly controlled series of prohormones and proteases. The unique, patented 'Immobilized Antigen' technology of this angiotensin II immunometric assay allows reliable detection of 1-2 pg/ml, or as little as 10% of the normal human plasma concentration.





#### ApoAl Monoclonal Antibody (Clone CC3821C4)

Lyophilized IgG<sub>1</sub> **Stability:** ≥1 year at 4°C

Summary: Antigen: human ApoAI amino acids 188-199 • Host: mouse, clone CC3821C4 • Cross Reactivity: (+) human ApoAI, (-) ApoB • Application(s): WB • ApoAI is a major protein component of HDL. It acts as an acceptor for sequential transfers of phospholipids and free cholesterol from peripheral tissues and transports cholesterol to the liver and other tissues for excretion and steroidogenesis.

#### **ApoAl Polyclonal Antibody**

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human ApoAI protein amino acids 188-199 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat ApoAI protein • Application(s): ICC and WB • ApoAI is a major protein component of HDL. It acts as an acceptor for sequential transfers of phospholipids and free cholesterol from peripheral tissues and transports cholesterol to the liver and other tissues for excretion and steroidogenesis.

• Also Available: ApoAI Blocking Peptide (13079)

#### ApoAl Western Ready Control

**Stability:** ≥6 months at -80°C

Summary: Source: human recombinant protein • Application(s): positive control

1 ea

#### 20-carboxy Arachidonic Acid

10007912

10009751

[79551-84-1] 20-COOH-AA MF:  $C_{20}H_{30}O_4$  FW: 334.5 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A major metabolite of 20-HETE produced in renal tubular epithelial, endothelial, and microvascular smooth muscle cells; inhibits ion transport in the kidneys and has vasodilatory activity

25 µg 50 µg

100 µg 500 µg

<sup>+</sup> SPI-BIO Assays are available through Cayman Chemical only within

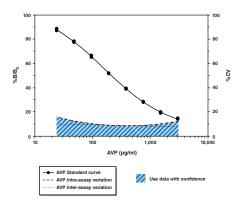
583951

ADH, Antidiuretic Hormone, Argipressin, AVP, Vasopressin

**Stability:** ≥1 year at -20°C **Sensitivity:** 50% B/B<sub>0</sub>: ~220-380 pg/ml • 80% B/B<sub>0</sub>: ~45-60 pg/ml

Summary: AVP is a nine amino acid peptide hormone that plays a primary role in the regulation of renal water excretion and a secondary role in the regulation of cardiovascular function in mammals. Cayman's AVP EIA is a competitive assay that can be used for the measurement of AVP from plasma and serum.

96 strip/solid wells 480 strip/solid wells



#### **Aspalatone**

[147249-33-0] Acetylsalicylic Acid Matol ester

MF:  $C_{15}H_{12}O_6$  FW: 288.3 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An anti-platelet aggregator (IC<sub>50</sub> = 180 µM, in vitro) that prolongs bleeding time significantly in a rodent model of thromboembolism; at 24 mg/kg, generates antioxidant and neuroprotective effects against kainic acidinduced epilepsy in rat hippocampus

5 mg 10 mg 50 mg 100 ma

**Aspirin** 

70260

[50-78-2] Acetylsalicylic Acid

**MF:** C<sub>0</sub>H<sub>8</sub>O<sub>4</sub> **FW:** 180.2 **Purity:** ≥99%

A crystalline solid **Stability:** ≥2 years at room temperature

Summary: A non-selective, irreversible COX inhibitor with IC<sub>50</sub> values of 0.75 and 1.25 mM for ovine COX-1 and COX-2, respectively

5 g 25 g 50 g 100 g

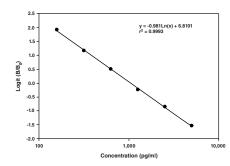


#### Aspirin™ Effect-Detection Kit

**Stability:** ≥1 year at 4°C

Summary: Cayman's Aspirin<sup>TM</sup> Effect-Detection assay is a 510K, clinically-approved diagnostic kit for the measurement of 11-dehydro TXB<sub>2</sub>. It is intended to help physicians assess the effectiveness of their patients' Aspirin™ regime, and to help identify the high-risk group who are inadequately controlled on an 80 mg aspirin dose. This assay can be completed in 3 hours and utilizes urine as the sample matrix. The assay exhibits intra-assay %CV values of <11% with sensitivity sufficient to detect the lower levels of 11-dehydro TXB<sub>2</sub> in patients that respond well to Aspirin™.

1 ea



#### Atherosclerosis Product Pack

10005292

10010153

**Stability:** ≥6 months at -20°C

Summary: Contains ACAT-1 polyclonal antibody and blocking peptide, cholesteryl linoleate hydroperoxides, POV-PC, and PGPC

1 ea

#### Atriopeptin (rat) EIA Kit+

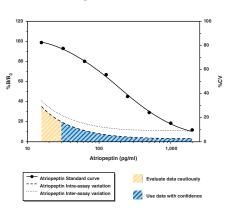
589401

**Stability:** ≥6 months at -20°C

**Sensitivity:** 50% B/B<sub>0</sub>: 190 pg/ml • 80% B/B<sub>0</sub>: 60 pg/ml

Summary: Atriopeptin is a 28 amino acid peptide synthesized primarily in cardiac atria. This peptide hormone acts in opposition to angiotensin II in regulating renal, hemodynamic, and endocrine function. Atriopeptin is released in response to the increased pressure and mechanical stretch of the right atrium due to blood volume overload. Atriopeptin then acts at the nephron to increase salt and water excretion, lowering blood volume and blood pressure.

96 wells



#### **AUDA** 10007927

[479413-70-2]

MF:  $C_{23}H_{40}N_2O_3$  FW: 392.6 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An inhibitor of sEH exhibiting IC<sub>50</sub> values of 18 and 69 nM for the murine and human enzymes, respectively

5 mg 10 mg 50 mg 100 mg

Azelaovl PAF

MF:  $C_{33}H_{66}NO_9P$  FW: 651.9 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A PAF analog isolated and purified from oxLDL; acts as a potent PPARy agonist

500 µg 1 mg 5 mg 10 mg

#### BAY-41-8543 10011131

[256498-66-5]

MF:  $C_{21}H_{21}FN_8O$  FW: 420.4 Purity:  $\geq$ 98%

A crystalline solid Stability: ≥2 years at -20°C

Summary: A stimulator of sGC, increasing the activity of recombinant sGC dosedependently, from 0.1 nM to 100 µM, up to 92-fold; in vitro, relaxes vessels and inhibits platelet aggregation at nM concentrations; in vivo, decreases blood pressure dose-dependently, prolongs bleeding time, and reduces thrombosis

#### 1-Benzylimidazole

[4238-71-5]

**MF:**  $C_{10}H_{10}N_2$  **FW:** 158.2 **Purity:**  $\geq$ 99%

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A selective inhibitor of TX synthase 1 g 5 g 25 g

100 g

#### Beraprost (sodium salt)

[88475-69-8] ML 1129, Procylin, TRK 100

**MF:**  $C_{24}H_{29}O_5$  • Na **FW:** 420.5 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An analog of prostacyclin in which the unstable enol-ether moiety has been replaced by a benzofuran ether function; exhibits a plasma half-life of several hours; improves survival and pulmonary hemodynamics in patients with primary pulmonary hypertension

1 mg 5 mg 10 mg 50 ma

#### Berberine 10006427

[633-65-8] BBR, Umbellatine

**MF:**  $C_{20}H_{18}ClNO_4$  **FW:** 371.8 **Purity:**  $\geq$ 95%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An alkaloid natural product that reduces total cholesterol, LDL cholesterol, and triglycerides in humans and hamsters; enhances LDL-receptor

protein and mRNA levels in hepatocytes

1 g 5 g 10 g 25 g

Bezafibrate

[41859-67-0] Benzofibrate, BM 15075, Bezalip, Bezatrol, Difaterol MF:  $C_{19}H_{20}CINO_4$  FW: 361.8 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A well-established pan PPAR agonist; activates human PPARα, PPARδ, and PPARy with EC50 values of 50, 20, and 60 µM, respectively; helps lower LDL cholesterol and triglycerides while raising HDL cholesterol levels

500 mg 1 g 5 g 10 g

#### BIBB 515

60924

70510

18230

10010517

10155

60928

[1566.35-05-1]

**MF:**  $C_{22}H_{21}ClN_2O_2$  **FW:** 380.9 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A selective and potent inhibitor of 2,3-oxidosqualene cyclase (OSC) in vivo with an ED<sub>50</sub> value of 0.2-0.5 and 0.36-33.3 mg/kg in rats and mice, respectively; reduced total cholesterol by 19% in normolipemic hamsters at a dose of 55 mg/kg/day for 11 days and 25% in hyperlipemic hamsters at a dose of 148 mg/kg/day for 25 days

5 mg 10 ma 25 ma

#### BM 567 [284464-77-3]

MF:  $C_{18}H_{28}N_4O_5S$  FW: 412.5 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A dual acting antithrombogenic agent; inhibitor of TXA<sub>2</sub> synthase (IC<sub>50</sub> = 12 nM); acts as a TP receptor antagonist ( $IC_{50} = 1.1 \text{ nM}$ )

NOTE: Sold under license from Université of Liège

#### **Butanoyl PAF**

1 mg

5 mg

10 mg

50 mg

**MF:** C<sub>28</sub>H<sub>58</sub>NO<sub>7</sub>P **FW:** 551.7 **Purity:** ≥98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A PAF analog isolated from oxLDL; retains at least 10% of the agonist potency of PAF itself but is present in oxLDL in amounts more than 100 times greater than PAF

1 mg 5 mg 10 mg 50 mg

#### **Butenoyl PAF** 60929 **MF:** C<sub>28</sub>H<sub>56</sub>NO<sub>7</sub>P **FW:** 549.7 **Purity:** ≥98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A PAF analog isolated from oxLDL; retains at least 10% of the agonist potency of PAF itself but is present in oxLDL in amounts more than 100 times

1 mg 5 mg 10 mg 50 mg

+ SPI-BIO Assays are available through Cayman Chemical only within North or South America and Asia: elsewhere contact SPI-BIC

CRP

10009893

C-Reactive Protein (human) EIA Kit

10011236

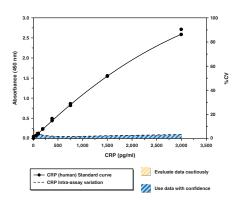
**Stability:** ≥6 months at 4°C

Limit of Detection: 46.9 pg/ml

Summary: CRP is a 224 amino acid protein that is synthesized primarily by hepatocytes, and to a lesser extent adipocytes. CRP plasma levels increase ~1,000fold in response to acute and chronic inflammatory conditions, making it a useful gauge of inflammation in a wide range of physiological and pathological conditions. Cayman's CRP (human) EIA is a sensitive immunometric assay which can be used to measure CRP in plasma without prior sample purification.

caymanchem.com

96 wells



#### Carbaprostacyclin

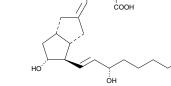
[69552-46-1] Carbacyclin, cPGI

**MF:**  $C_{21}H_{34}O_4$  **FW:** 350.5 **Purity:** ≥99%

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A stable analog of PGI2 with reduced potency; inhibits ADP-induced platelet aggregation in human PRP (ED<sub>50</sub> = 47 nM)

500 µa 1 mg 5 mg 10 mg



• Also Available: Carbaprostacyclin-biotin (18213)

Carbaprostacyclin methyl ester (9000183)

5-cis Carbaprostacyclin (18211)

#### 13,14-dehydro-15-cyclohexyl

Carbaprostacyclin [145375-81-1]

**MF:**  $C_{21}H_{30}O_4$  **FW:** 346.5 **Purity:** ≥98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A chemically stable analog of PGI<sub>2</sub>; inhibits the ADP-induced aggregation of human platelets with an ED<sub>50</sub> value of about 40 nM in PRP and 77 nM in washed platelets

50 µg 100 µg 500 µg

1 mg

#### Carbocyclic Thromboxane A<sub>2</sub>

[74034-56-3] Carbocyclic TXA<sub>2</sub>, CTA<sub>2</sub>

MF:  $C_{22}H_{36}O_3$  FW: 348.5 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

**Summary:** A stable analog of TXA<sub>2</sub> with potent coronary vasoconstrictor activity; inhibits arachidonic acid-induced aggregation (IC<sub>50</sub> = 4-5 μM); inhibits TXB<sub>2</sub> synthesis in rabbit platelets

100 µg 500 µg 1 mg 5 mg

#### Cardiogenol C (hydrochloride)

13187

10005186

19010

[671225-39-1]

**MF:**  $C_{13}H_{16}N_4O_2 \bullet HCl$  **FW:** 296.8 **Purity:**  $\geq$ 97%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: Induces the differentiation of MHC-positive cardiomyocytes from embryonic stem cells with an EC<sub>50</sub> value of 0.1 μM

1 mg 5 mg 10 mg 50 mg

CAY10441

18210

18212

[221529-58-4] RO1138452

MF:  $C_{10}H_{22}N_2O$  FW: 309.4 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A potent IP receptor antagonist; inhibits the binding of tritiated iloprost to rodent neuroblastoma cells with a K<sub>i</sub> value of about 1.5 nM

1 mg 5 mg 10 mg 25 mg

#### CAY10449 10005913

**MF:**  $C_{19}H_{21}N_3O_2$  **FW:** 323.4 **Purity:**  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A potent IP receptor antagonist; inhibits the binding of tritiated iloprost to rodent neuroblastoma cells with a K<sub>i</sub> value of about 3 nM

1 mg 5 mg 10 ma 25 mg

#### CAY10485

[615264-62-5] 3.4-dihydroxy Hydrocinnamic Acid (L-Aspartic Acid dibenzyl ester) amide MF:  $C_{27}H_{27}NO_7$  FW: 477.1 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An inhibitor of human ACAT-1 and ACAT-2 with IC<sub>50</sub> values of 95 and 81 µM, respectively; inhibits copper-mediated oxidation of LDL by 91% at a concentration of 2 µM

5 mg 10 mg 50 ma 100 ma

+ SPI-BIO Assays are available through Cayman Chemical only within North & South America and Asia; elsewhere contact SPI-BIO.

CAY10486 10006452

[615264-52-3] 4-Hydroxycinnamic Acid (L-phenylalanine methyl ester) amide MF:  $C_{19}H_{19}NO_4$  FW: 325.4 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An inhibitor of human ACAT-1 and ACAT-2 with an IC50 value of approximately 60 µM for both enzymes; also inhibits copper-mediated oxidation of LDL by about 28% at a concentration of 3 µM

5 mg 10 mg 50 mg 500 mg

#### CAY10487

10006480

[778624-05-8] 3,4-Dihydrocinnamic Acid (L-alanine methyl ester) amide MF:  $C_{13}H_{15}NO_5$  FW: 265.3 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An inhibitor of fatty streak lesion formation; also inhibits coppermediated oxidation of LDL by about 75% at a concentration of 2  $\mu M$ 

10 mg 50 mg 500 mg

CAY10499 10007875

[359714-55-9]

MF:  $C_{18}H_{17}N_3O_5$  FW: 355.3 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent inhibitor of human hormone sensitive lipase ( $IC_{50} = 90 \text{ nM}$ )

5 mg 10 mg 25 mg

#### CAY10514 10009017

[868526-38-9] Methyl-8-hydroxy-8-(2-pentyl-oxyphenyl)-oct-5-ynoate MF:  $C_{20}H_{28}O_4$  FW: 332.4 Purity:  $\geq 98\%$ 

A solution in methyl acetate **Stability:** ≥1 year at -20°C

Summary: An aromatic analog of 8(S)-HETE; acts as a dual agonist of PPARα and PPARγ with EC<sub>50</sub> values of 0.17 and 0.64 μM, respectively

1 mg .COOCH<sub>o</sub> 5 mg 10 mg 50 mg

#### CD36 Monoclonal Antibody (Clone JC63.1) 188150

GPIIIb, GPIV, Hexarelin Receptor, oxLDL Receptor, Thrombospondin Receptor Purified mouse anti-CD36 IgA **Stability:** ≥1 year at -20°C

Summary: Antigen: adenovirus expressing full-length recombinant murine CD36 • Host: CD36 null mouse, clone JC63.1 • Isotype: IgA • Cross Reactivity: (+) murine, rat, and human CD36 • Application(s): FC and functional blocking • CD36 is a type-B scavenger receptor that is necessary for the formation of foam cells in atherosclerotic lesions.

100 µg 500 µg CD36 Monoclonal Antibody (Clone JC63.1) (azide free)

GPIIIb, GPIV, Hexarelin Receptor, oxLDL receptor, Thrombospondin Receptor IgA **Stability:** ≥1 year at 4°C

Summary: Antigen: recombinant adenovirus expressing full-length mouse CD36 • Host: CD36 null mouse, clone JC63.1 • Cross Reactivity: (+) human, murine, and rat CD36 • Application(s): FC, functional blocking, and ICC • CD36 is a type-B scavenger receptor that is necessary for the formation of foam cells and thereby atherosclerotic lesions.

100 µg 500 µg

#### CD36 Monoclonal FITC Antibody (Clone JC63.1)

10009870

GPIIIb, GPIV, Hexarelin Receptor, oxLDL Receptor, Thrombospondin Receptor IgA-FITC **Stability:** ≥1 year at -20°C

Summary: Antigen: recombinant adenovirus expressing full-length mouse CD36 • Host: CD36 null mouse, clone JC63.1 • Cross Reactivity: (+) murine, rat, and human CD36 • Application(s): FC and ICC • CD36 is a type-B scavenger receptor that is necessary for the formation of foam cells and thereby atherosclerotic lesions.

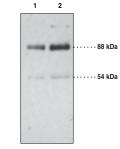
#### CD36 Polyclonal Antibody

100011

10007171

GPIIIb, GPIV, Hexarelin Receptor, oxLDL Receptor, Thrombospondin Receptor Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human CD36 amino acids 99-114 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat CD36 • Application(s): WB • CD36 is a type-B scavenger receptor that is necessary for the formation of foam cells in atherosclerotic lesions.



Lane 1: Human platelet lysate (15 µg)
Lane 2: Human platelet lysate (30 µg)

• Also Available: CD36 Blocking Peptide (300011)

#### Cetaben

[55986-43-1] Hexadecylamino-p-amino Benzoic Acid **MF:**  $C_{12}H_{20}NO_{2}$  **FW:** 361.6 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A unique, PPARa-independent peroxisome proliferator with hypolipidemic activity characterized by reduction in serum triglyceride and cholesterol concentrations in rats

10 mg 50 mg

# Epigenetics: The New Frontier for Vascular Diseases

In the world of personal healthcare, there aren't many things that seem as personally adjustable as cardiovascular health. There are 'heart smart' foods, suggesting that simple changes in diet can reap benefits for your heart. A regular schedule of aerobic exercise will strengthen your cardiovascular system. Adjust your sleep patterns, get a pet, find a spouse, eat less, or listen to classical music and you might enjoy a long life of vascular bliss. The ways

you can manipulate the system seem almost endless.

Naturally, there is the genetic side of vascular diseases. Genetic defects in such diverse parameters as lipid metabolism, arterial structure, endothelial and smooth muscle function, platelet adhesiveness, and fibrinolysis can clearly contribute to the development of atherosclerosis, for example. The phenotypic consequences of such defects may range from subtle to pronounced. Superficially, genetic problems seem so much more refractory to treatment than the personal, everyday lifestyle options.

Now, everything is about to get a lot more complicated as well as very interesting. The ways that all those adjustable, personal factors impact gene expression in prolonged, meaningful ways is part of the field of epigenetics. Technically, 'epigenetics' refers to heritable changes in gene expression that do not involve changes in DNA sequence or copy number. The cell biologist uses epigenetic events to help explain how pluripotent stem cells differentiate into diverse cell types, while developmental biologists see epigenetic processes at work in the morphogenesis of an embryo into discrete organs. Similarly, epigenetic signaling that is put in place in response to prenatal or early life events can contribute to disease development much later in life.

#### **Epigenetics: The Basics**

In molecular terms, epigenetic changes commonly involve adding or removing various "marks" on chromatin. The basic structural unit of chromatin is the nucleosome, which consists of a short stretch of DNA wrapped twice around a protein octamer composed of two sets of the core histones, H2A, H2B, H3, and H4 (Figure 1). The C-termini of the histones are highly conserved and form the protein core of each nucleosome, while the N-terminal tails project from the nucleosome. These tails can interact with DNA or with other proteins. For example, each tail is rich in regularly spaced, positively-charged residues (lysine (K), arginine (R)),



**Figure 1. Histone tails.** A.) Positioning of the histone tail relative to the C-terminal folded region. B.) Amino acid sequences of core histone N-terminal tails, indicating sites of phosphorylation (p), acetylation (a), ADP ribosylation (rib), and methylation (m).

which can associate with the anionic backbone of DNA. Acetylation or methylation of these residues alters their charge and size, affecting DNA binding. Phosphorylation, often on sites adjacent to basic residues, puts a large negatively-charged moiety on the histone. These types of chemical marks also lead to structural changes in the histone tail, providing binding sites for proteins or protein complexes, which in turn modulate further protein or histone changes. Histone marks can be reversed enzymatically, or they can persist for longer periods, extending through multiple cell divisions to maintain cell phenotype. A much more stable modification involves methylation of cytosine residues on DNA, most notably on 'CpG islands', genomic regions that are rich in CG pairs. Methylation of DNA is particularly stable and often results in the suppression of transcription of specific genes by physically interfering with the transcriptional machinery. Methylated DNA, like marked histones, provides binding sites for proteins and the establishment of protein complexes that can modify histones and alter gene expression. These are some of the basic ways that phenotype can be stably altered without changing DNA sequence or copy number.

#### Food For Thought

You are what you eat, but does your diet affect your children's health? About ten years ago, Bygren and colleagues studied a cohort of individuals born in 1905 in the Överkalix parish in northernmost Sweden, where annual harvests are heavily impacted by the weather. 1 County records provided birth and death dates for the 1905 cohort, their parents and their grandparents. Additional records indicated years of poor, moderate, or superior availability of crop food during the preceding century. Was the lifespan of the individuals born in 1905 affected by the nutritional experience of their predecessors? To focus this question, the authors hypothesized that, in order for famine or food surplus to have persistent effects, the dietary impact must occur in a sexually formative period. Interestingly, they found a significant correlation between food availability for paternal grandfathers when they were 9-12 years old and the survival of their grandchildren. Perhaps more remarkably, grandchild lifespan shortened if there was an excess of food for the paternal grandfather and increased if the grandfather experienced famine during this critical developmental stage. The effect was not trivial: the difference in survival for grandchildren averaged 32 years!

A follow-up study by the same group expanded the study to include cohorts from 1890, 1905, and 1920 and examined cause of death data.<sup>2</sup> They reported that a father experiencing famine during the critical 9-12 years of age passed protection against cardiovascular disease to his children. To restate: famine for the father meant less cardiovascular disease for his children. No significant correlations were found for mothers, grandmothers, or grandparents and offspring cardiovascular disease. These correlative studies anticipated a decade of studies into the now popular concept that the diet of the mother might reprogram the genes of the fetus, without mutation, leading to persistent effects in health and development of the child.<sup>3</sup>

#### Case in Point: Epigenetics in Hypertension

Aldosterone is a corticosteroid hormone that, through mineralocorticoid receptors, alters the uptake and secretion of salts by collecting ducts of the kidneys, thus altering water uptake, blood volume, and blood pressure. Excess aldosterone contributes to arterial hypertension, congestive heart failure, chronic kidney disease, coronary artery disease, and stroke.<sup>4</sup> Clearly, the mechanism of action of aldosterone at the kidney collecting duct must be important.

One enzyme serves to hypermethylate the lysine residue 79 on histone 3 (H3K79). In yeast, ablation of this methyltransferase causes disruption of telomere silencing, leading to the name DOT1. A mouse DOT1 homolog,

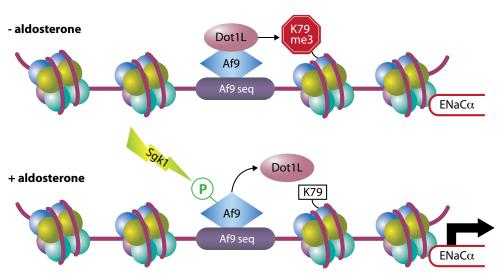


Figure 2. Aldosterone promotes the phosphorylation of Af9 by Sgk1, dissociating DOT1L from Af9, allowing demethylation of H3K79 and the expression of ENaCa.

DOT1L, also hypermethylates H3K79 in mice, but this serves to repress the expression of a subset of genes. DOT1L binds to AF9, a DNA-binding protein which positions DOT1L only on those nucleosomes whose associated DNA has a sequence that is recognized by AF9. This confers selectivity for which nucleosomes are targeted for hypermethylation of H3K79 and which genes are consequently repressed. One such gene encodes a sodium channel subunit, epithelial Na $^*$  channel subunit  $\alpha$  (ENaC $\alpha$ ).

Aldosterone alters the expression of ENaC $\alpha$  by modulating the DOT1L/AF9 repressive complex (Figure 2). Aldosterone induces the expression of serum- and glucocorticoid-induced kinase 1 (Sgk1), which phosphorylates AF9 and disrupts the complex.<sup>5</sup> Aldosterone also inhibits the expression of both DOT1L and AF9.<sup>6,7</sup> These steps are necessary and sufficient to increase ENaC $\alpha$  expression in the presence of the hormone-receptor complex.

It is worth noting that this pathway involves a short term and reversible mechanism involving histone modification. Such transient changes are the hallmark of the 'signaling model' of epigenetic modifications. These contrast with other pathways where histone modifications are inherited through cell division, the so-called 'histone code hypothesis' of epigenetic signaling.

#### **Classical Epigenetic Studies**

The Dutch Famine Birth Cohort Studies relate to babies born around a 5-month period of extreme food shortage during the winter of 1944-1945. Numerous characteristics of mothers and offspring were obtained around birth and the offspring have been followed since. In one study, exposure to famine during gestation led to a higher cumulative incidence of coronary artery disease, as well as an earlier age at onset, than was found for those not exposed to famine during gestation. § Moreover, these effects were found only if exposure to famine occurred early in gestation. Such correlative studies leave us to wonder: what could have happened during those first few months of development that became evident only years later?



Figure 3. Genetically identical week 15 littermates representing coat colors ranging from agouti (left) to pseudoagouti (right); Note differences in size

A series of recent studies have come closer to demonstrating epigenetics in action. While obviously the diet of a pregnant mother can affect fetal development, these studies indicate that dietary supplements, taken by the mother, may mark the genome of the fetus and affect adult health. In the viable yellow agouti (Avy) mouse, expression of the agouti gene leads to a switch from brown to yellow coat color (Figure 3). Moreover, adult agouti mice suffer from obesity, hyperlipidemia, and hypertension. The expression of the agouti gene is initiated from a cryptic promoter in a retrotransposon inserted in agouti pseudoexon 1A (PS1A).9 It's known that cytosine methylation on the transposable element prevents agouti gene expression, producing a brown-coated mouse that is referred to as 'pseudoagouti'. 10 Waterland and Jirtle demonstrated that supplementing normal chow given to pregnant mice with methyl donors (folic acid, betaine, vitamin B12, choline) produced an increase in methylation of CpG sites within PS1A in the offspring and shifted the coat color distribution toward the brown (pseudoagouti) phenotype.9 These methylation patterns were found in cells from diverse tissues, showing that Avy methylation is determined in the early embryo and maintained with high fidelity throughout development.

In another study, genistein, an isoflavonoid naturally found in soy, increased both PS1A methylation and frequency of pseudoagouti expression in offspring when added to the mother's diet.<sup>11</sup> Importantly, ectopic agouti expression is associated with adult-onset obesity, diabetes, and cancer.<sup>12</sup> Hypermethylation of PS1A, in mice from genistein-fed mothers, persisted into adulthood, decreasing agouti expression, and, remarkably, protecting mice from obesity.<sup>11</sup> Conversely, hypomethylation of PS1A, following exposure to the common chemical bisphenol A (BPA), increased agouti expression;<sup>13</sup> BPA is known to promote obesity, hyperlipidemia, and hypertension in mice. Taken together, these results demonstrate that maternal diet has in utero effects on the epigenome of the early embryo that can alter susceptibility to disease into adulthood.

#### Reference

- Bygren, L.O., Kaati, G., and Edvinsson, S. *Acta Biotheor.* **49**, 53-59 (2001). Kaati, G., Bygren, L.O., and Edvinsson, S. *Eur. J. Hum. Genet.* **10**, 682-688 (2002).
- Heerwagen, M.J., Miller, M.R., Barbour, L.A., et al. Am. J. Physiol. Regul. Integr. Comp. Physiol. 299(3), R711-R722 (2010).
- Treet wagen, M.S., Miller, M.R., Barbour, E.A., et al. Altr. S. Physiol. Negal. Integr. 1
   Calhoun, D.A. Circulation 114, 2572-2574 (2006).
- Zhang, W., Xia, X., Reisenauer, M.R., et al. J. Clin. Invest. 117(3), 773-783 (2007).
- Zhang, W., Xia, X., Reisenauer, M.R., et al. J. Biol. Chem. 281(26), 18059-18068 (2006).
- 7. Zhang, D., Yu, Z.-Y., Cruz, P., et al. Kidney Int. **75(3)**, 260-267 (2009).
- Painter, R.C., de Rooij, S.R., Bossuyt, P.M., et al. Am. J. Clin. Nutr. **84**, 322-327 (2006). Waterland, R.A. and Jirtle, R.L. *Mol. Cell. Biol.* **23(15)**, 5293-5300 (2003).
- Waterland, R.A. and Jirtle, R.L. Mol. Cell. Biol. 23(15), 5293-5300 (2003).
   Morgan, H.D., Sutherland, H.G.E., Martin, D.I.K., et al. Nat. Genet 23. 314-318 (1999).
- Morgan, A.D., Sutherland, A.G.E., Martin, D.I.N., et al. Nat. Genet. 25, 514-516 (1999).
   Dolinov D.C. Weidman, I.R. Waterland R.A. et al. Environ. Health Perspect. 114(4), 567-572 (2006).
- 12. Yen, T.T., Gill, A.M., Frigeri, L.G., et al. FASEB J. **8,** 479-488 (1994).
- 13. Dolinoy, D.C., Huang, D., and Jirtle, R.L. *Proc. Natl. Acad. Sci. USA* **104(32)**, 13056-13061 (2007).

CG 4305

18216

For a full product listing, please visit www.caymanchem.com

**MF:**  $C_{25}H_{32}O_4$  **FW:** 396.5 **Purity:** ≥95%

A solution in methyl accetate **Stability:** ≥1 year at -20°C

Summary: A stable carbacyclic PGI2 analog that inhibits ADP-induced platelet aggregation 25% at a concentration of 50 nM; 10 mg/kg (oral) or 1 mg/kg (intraduodenal) prevents thrombotic arterial occlusion

100 µg 500 µg 1 mg 5 mg

#### Chenodeoxycholic Acid

10011286

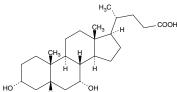
10012447

[474-25-9] Anthropodeoxycholic Acid, CDCA, Fluibil, Hekbilin, Kebilis, Ulmenide **MF:**  $C_{24}H_{40}O_4$  **FW:** 392.6 **Purity:**  $\geq$ 95%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A bile acid and farnesoid X receptor (FXR) ligand (EC<sub>50</sub> = 13-34  $\mu$ M) that is a key regulator of cholesterol homeostasis; exhibits toxicity that is linked to increased glutathione and increased oxidative stress; excess CDCA contributes to liver and intestinal cancers

1 g 5 g 10 g 25 g



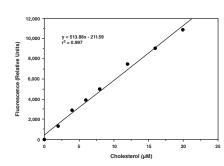
#### Cholesterol Assav Kit

10007640

**Stability:** ≥1 year at -20°C

Summary: Cholesterol, particularly in the form of LDLs, is well understood to be associated with increased risk of coronary heart disease. The measurement of cholesterol is one of the most common tests performed in the clinical laboratory setting. However, simple and easy assays for cholesterol in the research lab have not been readily available. Cayman's Cholesterol Assay provides a simple fluorometric method for the sensitive quantitation of total cholesterol in plasma or serum.

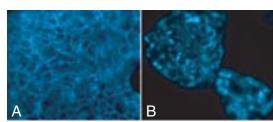
96 wells 480 wells



#### Cholesterol Cell-Based **Detection Assay Kit**

**Stability:** ≥6 months at -20°C

Summary: The mechanism for the movement of cholesterol from intracellular sites to their ultimate cellular destination is an unresolved question of fundamental importance to cell biology and medicine. Cayman's Cholesterol Cell-based Detection Assay provides a simple fluorometric method to study mechanisms and biological factors that regulate cholesterol metabolism or movement within cells.



Accumulation of cholesterol inside HepG2 cells in response to 1.25 μM U18666A. HepG2 cells were treated with DMSO (vehicle) or 1.25  $\mu M$ U-18666A for 48 hours. Panel A: Cells treated with DMSO alone demonstrate that majority of cholesterol is localized on the plasma membrane. Panel B: U-18666A treatment for 48 hours induces intracellular accumulation of cholesterol droplets, indicating blockage of intracellular cholesterol transport.

#### 3-dodecanovl-NBD Cholesterol

13220

10009779

3-C12-NBD Cholesterol

MF:  $C_{45}H_{70}N_4O_5$  FW: 747.1 Purity:  $\geq 98\%$ 

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A fluorescently-tagged cholesterol with the hydrophilic NBD fluorophore attached to the hydrophilic end of cholesterol, separated by a 12-carbon spacer; allows the cholesterol to properly orient in membrane bi-layers while the fluorescent tag is presented outside of the bi-layer

mg 5 mg

#### 3-hexanoyl-NBD Cholesterol

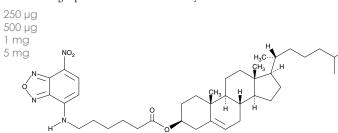
13221

3-C6-NBD Cholesterol

MF: C<sub>20</sub>H<sub>50</sub>N<sub>4</sub>O<sub>5</sub> FW: 662.9 Purity: ≥98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A fluorescently-tagged cholesterol with the hydrophilic NBD fluorophore attached to carbon 3, at the hydrophilic end of cholesterol, separated by a 6-carbon spacer; allows the cholesterol to properly orient in membrane bi-layers while the fluorescent tag is presented outside of the bi-layer



#### 5α-hydroxy-6-keto Cholesterol

[13027-33-3] Cholestane-6-0x0-3β,5α-diol, 6-Ox0-3,5-diol

**MF:** C<sub>27</sub>H<sub>46</sub>O<sub>3</sub> **FW:** 418.7 **Purity:** ≥98%

5 mg

10 mg

50 mg

100 µg

500 µg

1 mg

5 mg

Cicaprost

certain cell types

500 µg

1 mg

5 mg

10 mg

[94079-80-8] ZK 96480

MF:  $C_{22}H_{30}O_5$  FW: 374.5 Purity:  $\geq$ 98%

A solution in methyl acetate **Stability:** ≥1 year at -20°C

Summary: A prostacyclin analog that is orally active with prolonged availability in

vivo, having a terminal half-life in plasma of one hour; inhibits the pro-inflammatory

actions of certain leukocytes, suppresses cardiac fibrosis, and blocks mitogenesis of

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A major metabolite of cholesterol formed during exposure of lung epithelial cells to ozone; potent inhibitor of cholesterol synthesis in human bronchial epithelial cells (IC<sub>50</sub> = 350 nM); exhibits significant cytotoxicity in the low  $\mu$ M range

Summary: A product derived from the autoxidation of cholesteryl linoleate

Cholesteryl Linoleate Hydroperoxides

A solution in ethanol **Stability:** ≥6 months at -80°C

MF: C<sub>45</sub>H<sub>76</sub>O<sub>4</sub> FW: 681.1 Purity: ≥98% hydroperoxide content

containing a mixture of racemic 9- and 13-HpODE cholesteryl esters

10007601

48001

16831

5 mg

10 mg

50 mg

Clofibrate [637-07-0]

**MF:** C<sub>12</sub>H<sub>15</sub>ClO<sub>3</sub> **FW:** 242.7 **Purity:** ≥98% A colorless liquid **Stability:** ≥1 year at -20°C

Ciprostene (calcium salt)

**MF:**  $[C_{22}H_{36}O_4]_2 \cdot Ca^{2+}$  **FW:** 769.1 **Purity:**  $\ge 98\%$ 

A crystalline solid **Stability:** ≥6 months at -20°C

[81703-55-1] U-61431F

Summary: A PPARα agonist used clinically to treat dyslipidemia and cardiovascular disease; exhibits EC<sub>50</sub> values of 50 and 55 μM for murine and human PPARα,

Summary: A 9β-methyl analog of carbaprostacyclin and a stable analog of PGI<sub>2</sub>;

exhibits biological activity similar to PGI<sub>2</sub>, but is 30-fold less potent

500 µl 1 ml 5 ml 10 ml

#### (±)-Clopidogrel (hydrochloride)

13657

9000368

10005745

**MF:** C<sub>16</sub>H<sub>16</sub>NO<sub>2</sub>S • HCl **FW:** 358.3 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An antithrombic compound whose active metabolite is a selective, irreversible antagonist of the platelet purinergic P2Y receptor (IC<sub>50</sub> = 100 nM); inhibits ADP-induced platelet aggregation ex vivo; shown to be beneficial in the prevention of vascular ischemic events for patients without deficiencies in CYP2C19related metabolism

10 mg 50 mg

#### 3-hydroxy-3-methylglutaryl-

Coenzyme A-d<sub>3</sub> (ammonium salt)

MF:  $C_{27}H_{38}D_3N_7O_{20}P_3S \cdot 4NH_4$  FW: 983.8 Chemical Purity: ≥96%

**Deuterium Incorporation:** ≤1% d<sub>o</sub>

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: Intended for use as an internal standard for the quantification of HMG-CoA by GC- or LC-MS

100 µg 500 µg 1 mg 5 mg

#### For Laboratory Use Only. Not for human or veterinary diagnostic or therapeutic use. Prices are US retail list price as of 05-01-2011. For current US pricing

#### COX-1 Monoclonal Antibody (Clone CX111) 160110

Cyclooxygenase 1, PGHS-1, PGH Synthase 1

Lyophilized IgG **Stability:** ≥3 years at -20°C

**Summary:** Antigen: purified ovine COX-1 • Host: mouse, clone CX111 • Isotype: IgG<sub>28</sub> • Cross Reactivity: (+) ovine, bovine, human, murine, rat, and monkey COX-1, ovine COX-2 (50%), and human COX-2 (~5%); (-) murine COX-2 • Application(s): IHC and WB • Constitutively expressed COX-1 is responsible for the production of PGs essential for the normal function of many organs including stomach and kidney.

1 ea · 70 kDa

• Also Available: COX-1 Monoclonal FITC Antibody (Clone CX111) (160111)

COX-1 Monoclonal PE Antibody (Clone CX111) (160120)

Lane 1: Ovine COX-1 (microsomal fraction

COX-1 (murine) Polyclonal Antibody (160109)

COX-1 (ovine) Polyclonal Antiserum (160108)

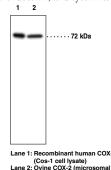
#### COX-2 Monoclonal Antibody (Clone CX229) 160112

Cyclooxygenase 2, PGHS-2, PGH Synthase 2

Lyophilized IgG Stability: ≥3 years at -20°C

Summary: Antigen: human COX-2 amino acids 580-599 • Host: mouse, clone CX229 • Isotype: IgG<sub>1</sub> • Cross Reactivity: (+) human, ovine, and monkey COX-2 (-) murine, rat, and rabbit COX-2 and COX-1 (all species) • Application(s): ICC, IHC, and WB • Expression of COX-2, the inducible form of COX, is regulated by growth factor tumors, tumor promoters, bacterial endotoxin, and cytokines.

1 ea



• Also Available: COX-2 (human) Blocking Peptide (360107)

COX-2 (human) Polyclonal Antibody (160107)

Goat Anti-COX-2 (human) Affinity-Purified

Polyclonal Antibody (100034)

COX-2 Monoclonal FITC Antibody (Clone CX229) (160113)

COX-2 (murine) Blocking Peptide (360106)

COX-2 (murine) Polyclonal Antibody (160106) COX-2 (murine) Polyclonal Antibody (160126)

COX-2 (murine) Polyclonal FITC Antibody (10010096)

COX-2 (murine) Polyclonal Antiserum (160116)

#### CUDA 10007923

[479413-68-8]

**MF:**  $C_{19}H_{36}N_2O_3$  **FW:** 340.5 **Purity:**  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** An inhibitor of sEH exhibiting IC<sub>50</sub> values of 11.1 and 112 nM for the murine and human enzymes, respectively

10 mg 25 mg 50 mg

#### N,N'-Dicyclohexylurea

1000497

[2387-23-7] DCU, NSC 17013

MF:  $C_{13}H_{24}N_2O$  FW: 224.3 Purity:  $\geq 98\%$ A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A selective sEH inhibitor with IC<sub>50</sub> values of 160 and 90 nM for recombinant human and murine sEH, respectively; at 10 µM, blocks the hydrolysis of cis- and trans-EETs and completely inhibits 14(15)-EET-induced PPAR activation

10 g 25 g 50 g 100 g		
25 g	0 11	
50 g	N N	H
100 g	Ĭ Ï	
	н 🖊	\

#### Diphenyl-1-pyrenylphosphine

62237

[110231-30-6] DPPP

**MF:** C<sub>28</sub>H<sub>19</sub>P **FW:** 386.4 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A probe that reacts stoichiometrically with hydroperoxides to yield the fluorescent molecule DPPP-O; also a fluorescent probe for the detection of LDL and cellular oxidation



#### 14,15-EE-8(Z)-E

10010486

[519038-93-8] 14,15-Epoxyeicosa-8(Z)-enoic Acid

MF:  $C_{20}H_{36}O_3$  FW: 324.5 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A structural analog of 14(15)-EET that demonstrates potent vasodilator agonist activity in bovine coronary arteries similar to that of 14(15)-EET

~	•	•	
25 µg			
60 µg			
00 110			<
00 µg			
600 µg			$\overline{}$
13			<b>*</b> 0 <b>*</b>

NOTE: relative stereochemistry shown in chemical structure

• Also Available: 14,15-EE-5(Z)-E (10004946)

#### Eicosapentaenoic Acid ethyl ester

10008884

[86227-47-6] EPA-EE, EPA ethyl ester

**MF:**  $C_{22}H_{34}O_2$  **FW:** 330.5 **Purity:**  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A stabilized ethyl ester form of this ω-3 C20:5 PUFA; dietary EPA-EE in rats increases fatty acid β-oxidation enzyme levels, down-regulates lipogenic genes, and decreases plasma cholesterol and triglyceride levels; blocks induced insulin resistance and corrects changes in adiponectin levels and TNF-α expression in rats fed a high-fat diet



• Also Available: Eicosapentaenoic Acid (90110)

Eicosapentaenoic Acid-d<sub>s</sub> (10005056)

Eicosapentaenoic Acid (peroxide free) (90110.1)

()-3 Fatty Acids				
Item No.	Product Name	Sizes		
90011	ω-3 Arachidonic Acid	1 mg • 5 mg • 10 mg • 25 mg		
90310	Docosahexaenoic Acid	50 mg • 100 mg • 250 mg • 500 mg		
90165	Docosapentaenoic Acid	1 mg • 5 mg • 10 mg • 100 mg		
90170	Docosatrienoic Acid	10 mg • 25 mg • 50 mg • 100 mg		
90110	Eicosapentaenoic Acid	50 mg • 100 mg • 250 mg • 500 mg		
90210	$\alpha$ -Linolenic Acid	50 mg • 100 mg • 250 mg • 1 g		
90320	Stearidonic Acid	1 mg • 5 mg • 25 mg		

For a full product listing, please visit www.caymanchem.com

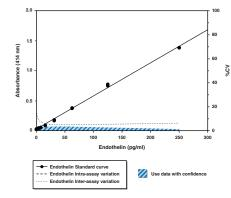
#### **Endothelin EIA Kit**

583151

**Stability:** ≥6 months at -20°C Limit of Detection: 7.8 pg/ml

Summary: The endothelin peptide family consists of three isoforms, ET-1 (corresponding to the initially isolated and most predominant isoform), ET-2, and ET-3. ET-1 is a 21 amino acid peptide and is one of the most potent vasoconstrictors currently known. ET-2 displays similar pharmacology to ET-1, whereas ET-3 is a weak vasoconstrictor but more potent inhibitor of platelet aggregation. Cayman's Endothelin Assay is an immunometric (i.e., sandwich) EIA that permits endothelin measurements within the range of 0-250 pg/ml, typically with a limit of detection of 1.5 pg/ml. Inter- and intra-assay CV's of less than 10% may be achieved at most concentrations. This assay offers sensitive and specific analysis of endothelin in serum, plasma, urine, or cell culture media.

96 wells 480 wells

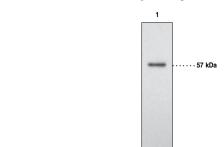


#### Endothelial Lipase Polyclonal Antibody

EDL, EL

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human endothelial lipase amino acids 19-32 • Host: rabbit • Cross Reactivity: (+) human, murine, rat, porcine, and ovine endothelial lipase • Application(s): IHC and WB • EL is a major genetic determinant for the concentration, structure, and metabolism of HDL, which protects against atherosclerosis.



Lane 1: HepG2 cell lysate (~30 µg)

• Also Available: Endothelial Lipase (human) Blocking Peptide (10004111)

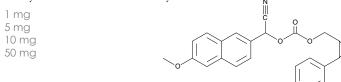
#### **Epoxy Fluor 7**

[863223-43-2]

**MF:** C<sub>23</sub>H<sub>19</sub>NO<sub>5</sub> **FW:** 389.4 **Purity:** ≥98%

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A sensitive fluorescent substrate for sEH that can be used to monitor the activity of both human and murine enzymes



#### Fenofibrate

10005368

70440

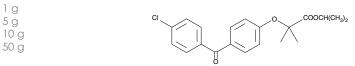
10008610

[49562-28-9]

MF:  $C_{20}H_{21}ClO_4$  FW: 360.8 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A PPARa agonist and hypolipidemic drug used clinically to treat dyslipidemia and cardiovascular disease; exhibits EC<sub>50</sub> values of 18 and 30 µM for murine and human PPARα, respectively



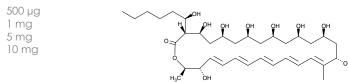
#### Filipin III

[480-49-9]

MF:  $C_{35}H_{58}O_{11}$  FW: 654.8 Purity:  $\geq 85\%$ 

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A fluorescent molecule that can be used to detect cholesterol by fluorescence and freeze-fracture electron microscopy



#### Fluvastatin (sodium salt)

10010337

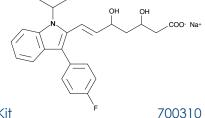
[93957-55-2]

MF:  $C_{24}H_{25}FNO_4$  • Na FW: 434.4 Purity: ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent, competitive inhibitor of HMG-CoA reductase (K<sub>i</sub> = 0.3 nM)

10 mg 25 mg 50 mg

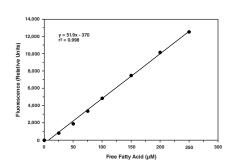


#### Free Fatty Acid Assay Kit

**Stability:** ≥6 months at -20°C

Summary: The measurement of FFA can be useful in determining metabolic status. Cayman's FFA Assay provides a simple, reproducible, and sensitive tool for measuring FFA's in plasma, serum, and urine. The FFA Assay utilizes a coupled enzymatic reaction that results in generation of the highly fluorescent product resorufin.

96 wells



#### Furegrelate (sodium salt)

[85666-17-7] U-63557A

MF:  $C_{15}H_{10}NO_3 \cdot Na$  FW: 275.2 Purity:  $\geq 99\%$ A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent inhibitor of TX synthase ( $IC_{50} = 15 \text{ nM}$ )

5 mg 10 mg 50 mg

GSK264220A

[685506-42-7]

MF:  $C_{17}H_{21}N_3O_4S$  FW: 363.4 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent inhibitor of endothelial lipase ( $IC_{50} = 16 \text{ nM}$ )

5 mg 10 mg 25 mg

#### Guanylate Cyclase $\alpha$ subunit (soluble) Polyclonal Antibody

sGC α subunit

Affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human guanylate cyclase α subunit amino acids 418-436 • Host: rabbit • Cross Reactivity: (+) murine, human, and bovine guanylate cyclase α subunit • Application(s): WB • Soluble guanylate cyclase is a heterodimeric enzyme, composed of  $\alpha$  and  $\beta$  subunits, that synthesizes cGMP from GTP. The enzyme is activated by the binding of nitric oxide or carbon monoxide to the heme group of the enzyme.

500 ul

• Also Available: Guanylate Cyclase α, subunit (soluble) Blocking Peptide (360895)

#### Guanylate Cyclase β<sub>1</sub> subunit (soluble) Polyclonal Antibody

160897

160895

sGC β, subunit

Affinity-purified IgG **Stability:** ≥1 year at -20°C

**Summary:** Antigen: rat soluble guanylate cyclase  $\beta_1$  subunit amino acids 188-207 • Host: rabbit • Cross Reactivity: (+) most mammalian species • Application(s): IHC and WB  $\bullet$  Soluble guanylate cyclase is a heterodimeric enzyme, composed of  $\alpha$  and  $\beta$ subunits, that synthesizes cGMP from GTP. The enzyme is activated by the binding of nitric oxide or carbon monoxide to the heme group of the enzyme.

500 ul

• Also Available: Guanylate Cyclase β<sub>1</sub> subunit (soluble) Blocking Peptide (360897)

#### (E)-Guggulsterone

10011296

[39025-24-6]

MF:  $C_{21}H_{28}O_2$  FW: 312.5 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A competitive antagonist of FXR (IC<sub>50</sub> = 15  $\mu$ M) that inhibits chenodeoxycholic acid (CDCA)-induced transactivation of FXR; lowers LDL cholesterol and triglyceride levels in rodents fed a high cholesterol diet

5 mg 10 mg 25 mg 50 mg

70540

#### (Z)-Guggulsterone

71800

[39025-23-5]

5 mg

10 mg

25 mg

MF:  $C_{21}H_{28}O_2$  FW: 312.5 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A competitive antagonist of FXR (IC<sub>50</sub> = 17  $\mu$ M ) that lowers LDL cholesterol and triglyceride levels in rodents fed a high cholesterol diet; demonstrates antitumor-promoting effects in human multiple myeloma and DU145 human prostate cancer cells

#### GW 0742 10006798

[317318-84-6]

**MF:**  $C_{21}H_{17}F_4NO_3S_2$  **FW:** 471.5 **Purity:**  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A selective PPAR $\delta$  agonist (EC<sub>50</sub> = 1.1 nM) that exhibits 1,000-fold selectivity over the other human PPAR subtypes

5 mg 10 mg 25 mg 50 mg

#### GW 7647 10008613

[265129-71-3]

**MF:**  $C_{29}H_{46}N_2O_3S$  **FW:** 502.8 **Purity:**  $\geq$ 98%

A crystalline solid Stability: ≥2 years at -20°C

**Summary:** A potent, selective PPARα agonist; activates human PPARα, PPARγ, and PPARδ with EC<sub>50</sub> values of 0.006, 1.1, and 6.2 μM, respectively

GW 9578 10011211

[247923-29-1]

**MF:**  $C_{26}H_{34}F_2N_2O_3S$  **FW:** 492.6 **Purity:**  $\geq$ 98%

A solution in methyl acetate **Stability:** ≥1 year at -20°C

**Summary:** A potent, selective PPARα agonist that activates the murine and human receptor with EC<sub>50</sub> values of 0.005 and 0.05 μM, respectively; 0.2 mg/kg given orally once daily for three days decreased serum total LDL cholesterol 40-60% in male Sprague-Dawely rats

#### GW 590735 10009880

[622402-22-6]

**MF:**  $C_{23}H_{21}F_3N_2O_4S$  **FW:** 478.5 **Purity:**  $\geq$ 98% A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent and selective PPAR $\alpha$  agonist (EC<sub>50</sub> = 4 nM) that is at least 500fold selective for PPARα over PPARγ and PPARδ

1 mg 5 mg 10 mg 25 mg

#### Hesperetin 10006084

[520-33-2]

MF:  $C_{16}H_{14}O_6$  FW: 302.3 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A citrus flavonoid that lowers plasma cholesterol; reduces the transcription of ACAT-2 mRNA in HepG2 cells and reduces ApoB protein synthesis (EC $_{50}$  ~ 50  $\mu$ M)

25 g 50 g 100 a 500 g 75780 HET0016

[339068-25-6] N-hydroxy-N'-(4-n-butyl-2-methylphenyl)Formamidine

**MF:**  $C_{12}H_{18}N_2O$  **FW:** 206.3 **Purity:**  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An inhibitor of 20-HETE formation in human renal microsomes with an IC<sub>50</sub> value of 8.9 nM, selectively inhibiting CYP4A and 4F isoforms

1 mg 5 mg 10 mg 100 ma

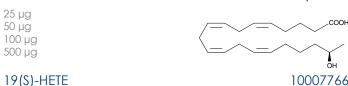
19(R)-HETE 10007767

[115461-39-7]

MF:  $C_{20}H_{32}O_3$  FW: 320.5 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A major CYP450 metabolite of arachidonic acid that is released from the kidney in response to angiotensin II; a potent vasodilator of renal preglomerular vessels; blocks 20-HETE-induced vasoconstriction of renal arterioles at 1 µM



[115461-40-0]

**MF:**  $C_{20}H_{32}O_3$  **FW:** 320.5 **Purity:**  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A major CYP450 metabolite of arachidonic acid that is released from the kidney in response to angiotensin II; a potent vasodilator of renal preglomerular

25 µg 50 µg 100 µg 500 µg

#### (±)9-HODE cholesteryl ester

38401

38601

[33783-76-5]

25 ua

50 µg

100 µg

250 µg

MF:  $C_{45}H_{76}O_3$  FW: 665.1 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A racemic monohydroxy fatty acid cholesteryl ester found in atherosclerotic lesions

• Also Available: 9(R)-HODE cholesteryl ester (38406) 9(S)-HODE cholesteryl ester (38411)

#### (±) 13-HODE cholesteryl ester

[167354-91-8]

**MF:**  $C_{45}H_{76}O_3$  **FW:** 665.1 **Purity:**  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A racemic monohydroxy fatty acid cholesteryl ester found in atherosclerotic lesions

25 µg 50 µg 100 µg 500 µg

Also Available: 13(R)-HODE cholesteryl ester (38606) 13(S)-HODE cholesteryl ester (38611) 20 Cayman Chemical caymanchem.com Cayman Chemical

#### Thomas G. Brock, Ph.

# You Say 'Thrombosis', I Say 'Thromboxane'



A blood clot that forms internally, in a blood vessel, looks different from the familiar scab that closes the laceration on your skin. Clots in the blood form lumps suggestive of curdling of the blood, if such a thing could happen. For this reason, it makes sense that such a clot is called a 'thrombus', after the Greek word thrombos, a lump or curd. Thrombosis, then, is simply the formation or presence of a blood clot in a blood vessel, which may be any vein or artery. An important early step in the formation of the thrombus is the aggregation of platelets at the blood vessel wall. One mediator that potently promotes platelet activation and aggregation is thromboxane A<sub>2</sub>. (TXA<sub>2</sub>). This article touches on the interrelationship of thromboxane and thrombi in certain types of thrombosis.

#### Thrombosis

The formation of the primary platelet plug in the damaged blood vessel is called primary hemostasis. The healthy endothelium inhibits platelet activation by secreting nitric oxide (NO) and prostaglandin I<sub>2</sub> (PGI<sub>2</sub>, or prostacyclin). However, disruption of the endothelial barrier through injury or disease exposes elements of the subendothelium to which platelets can adhere, including von Willebrand factor (vWf), glycoproteins, and extracellular matrix proteins (Figure 1). The binding of platelets is facilitated by fibrinogen. Adhesion of platelets to the vessel wall causes platelet activation, characterized by the release of pre-formed mediators from granules and the rapid synthesis of thromboxane A2 (TXA2) and platelet activating factor (PAF). These potently activate additional platelets and drive platelet aggregation, as well as vasoconstriction. With fibrinogen, aggregating platelets form a primary platelet plug, which is stabilized by the formation of fibrin through the coagulation cascade (a portion of secondary hemostasis). Continued coagulation produces a mesh of insoluble crosslinked fibrin, interspersed with plug remnants, that forms the core of the clot. Normally, clots are broken down by plasmin, which degrades the cross-linked fibrin. This depends on plasminogen activators, like tissue plasminogen activator (tPA), converting plasminogen to plasmin, which mediates fibrinolysis. Interestingly, many of the factors that are necessary for clot removal are produced by the healthy endothelium, so that the clot will not be removed until the initial damage to the endothelium is repaired.

The term 'hemostasis' literally refers to a balance, maintained in the blood, between clotting, anticoagulant, and fibrinolytic forces. Because the processes that contribute to this balance are complex, there are many genetic, acquired, and environmental factors that can tip the balance in favor of coagulation and cause the pathologic formation of thrombi. These can occur in arteries, veins, or cardiac chambers. Clots in arteries interfere with the delivery of oxygen to tissues, producing outcomes like myocardial infarction or ischemic stroke. Venous thrombosis commonly involves the activation of coagulation in vessels with reduced blood flow. The thrombus can interfere with blood flow at the site of formation, cause inflammation (phlebitis), or it can break free from the vessel wall and travel through the bloodstream as a thromboembolism, blocking a distant blood vessel. Death from deep vein thrombosis is usually due to massive pulmonary embolism, which causes some 300,000 deaths annually in the United States.

#### Thromboxane

Given the pivotal role of platelet plug formation in initiating clotting, current therapies focusing on the prevention of thrombosis seek to block platelet activation and aggregation. Clopidogrel blocks low affinity ADP receptors on platelets, preventing fibrinogen binding, short-circuiting the clotting process. While efficacious in a large percentage of individuals, there remain many for whom clopidogrel does not work. Moreover, the production of  $TXA_2$  can be increased in people with type II diabetes, <sup>1</sup> and clopidogrel does not solve this problem.

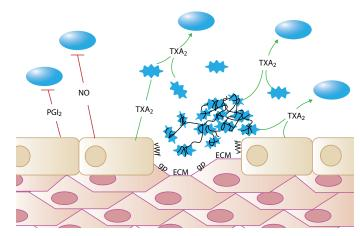


Figure 1. Healthy endothelium produces PGI<sub>2</sub> and NO, which suppresses platelet activation. Disruption of the endothelium reveals von Willebrand factor (WVf), glycoproteins (gp), and extracellular matrix (ECM), which induce platelets to change shape, aggregate, and release mediators, including TXA<sub>2</sub>. TXA<sub>2</sub> from activated platelets triggers additional platelet activation and aggregation.

TXA<sub>2</sub> is a lipid messenger derived by the enzymatic modification of arachidonic acid (AA), a polyunsaturated fatty acid that is stored in all cellular membranes, including those of platelets (Figure 2). Upon platelet activation, AA is released in abundance by cytosolic phospholipase A2 (cPLA2), an enzyme that specifically targets membrane phospholipids containing AA. Cyclooxygenase-1 (COX-1) and COX-2 isozymes catalyze the conversion of AA to prostaglandin (PG)H<sub>2</sub>, the initial step in the synthesis of TXA<sub>2</sub>, PGI<sub>2</sub>, and other primary prostaglandins. COX-1 is constitutively expressed in many cell types and, because of the stability of the protein, remains in platelets long after their formation from megakaryocytes. The expression of COX-2, on the other hand, is induced by inflammatory mediators and the enzyme typically has a relatively short half-life. Platelets typically lack COX-2. While most non-steroidal anti-inflammatory drugs (NSAIDs) reversibly inhibit both COX isoforms, aspirin inactivates them irreversibly. Coxibs selectively inhibit COX-2. Because platelets lack nuclei and cannot synthesize new COX-1, low-dose aspirin is sufficient to permanently inactivate COX-1 in platelets, preventing the formation of TXA2 that drives platelet activation and aggregation. This treatment is particularly effective in the prevention of arterial thrombi, which potentially contribute to heart attack and stroke. TXA2 synthase (TXAS) completes the synthesis of TXA2 from PGH<sub>2</sub>. TXA<sub>2</sub> has a half-life of about 30 seconds in plasma and is nonenzymatically hydrolyzed to TXB2, which is functionally inactive.

The focus above is largely on platelets as both the producers and responders of  $TXA_2$ . Importantly, other cells, including endothelial cells and leukocytes, can synthesize  $TXA_2$ . More importantly, neighboring cells can cooperate in the synthesis: if platelets do not use AA to make  $TXA_2$ , they can donate the AA to nearby cells, which can then use it to produce lipid mediators, including  $TXA_2$ ,  $PGI_2$ , and  $PGE_2$ . These products are then secreted, acting on platelets and adjacent cells. Such 'transcellular' synthesis may be particularly effective in the smaller vessels, where circulation slows and cells of different types are in close proximity.

#### Thromboxane Receptors

In general terms, most of the primary COX products, including  $TXA_2$ , activate G protein-coupled receptors (GPCR) at the cell surface. The binding pocket for GPCRs is typically deep in the membrane, defined by a few of the seven transmembrane domains of the receptor. Importantly, the binding pocket of each GPCR is highly selective for each ligand. As a result,

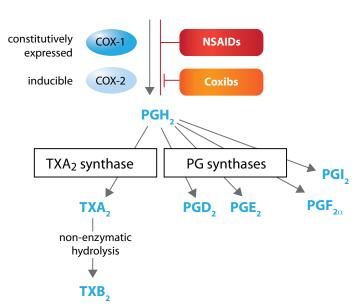
TXA2 binds one receptor, called the 'T prostanoid' (TP) receptor, which has two isoforms (TPα and TPβ). As with other GPCRs, agonist binding produces a change in receptor conformation, exposing intracellular sites involved in the interaction with the G protein. The G protein, consisting of  $\alpha$ ,  $\beta$ , and  $\gamma$  subunits, interacts with the receptor, leading to the release of guanosine diphosphate (GDP) from the Ga subunit and its replacement with guanosine triphosphate (GTP). The binding of GTP activates  $G\alpha$  so that it dissociates from the  $G\beta\gamma$  dimer and triggers a  $G\alpha$ -specific pathway. Both TP $\alpha$  and TP $\beta$  act primarily through G $\alpha$ q, which activates phospholipase CB, hydrolyzing phosphatidylinositol-4,5-bisphosphate (PIP<sub>2</sub>) to generate diacylglycerol (DAG) and inositol-1,4,5-trisphosphate (IP<sub>3</sub>). These mediators lead to a transient increase in intracellular calcium and activate certain isoforms of protein kinase C (PKC) and various calmodulindependent enzymes (e.g., kinases, NOS). Both TPα and TPβ can also act through Ga12-RhoA, activating Rho-associated protein kinase (ROCK), which is particularly important in regulating cytoskeletal components involved in contractility, mobility, and mitosis.

Both  $TP\alpha$  and  $TP\beta$  are derived from a single gene but are the product of unique mRNAs whose expressions are controlled by distinct promoters that are differentially regulated.<sup>2,3</sup> Structurally, the two receptor isoforms are identical except for the carboxy tails: fifteen residues at the terminus of  $TP\alpha$  are replaced with a much longer segment of 79 amino acids in  $TP\beta$  (Figure 3). These variants offer unique sites for phosphorylation and desensition of each isoform. For example, elevation of cytoplasmic cAMP levels activates protein kinase A (PKA), which phosphorylates  $Ser^{329}$  and  $Ser^{331}$  on  $TP\alpha$ , leading to desensitization of the receptor.<sup>4,5</sup> Different amino acids are present at these sites on  $TP\beta$ , so this receptor isoform is not affected by changes in cellular cAMP levels. Similarly,  $Ser^{331}$  is targeted by protein kinase G (PKG), which is activated by cGMP, an effector of NO, only on  $TP\alpha$ .<sup>4</sup> Interestingly, both isoforms are desensitized by PKC, although different residues are phosphorylated ( $Thr^{337}$  on  $TP\alpha$ ,  $Thr^{399}$  on  $TP\beta$ ).<sup>5</sup> Much more information on the TP isoforms is available.<sup>6</sup>

#### **Modulating Thromboxane Signaling**

The central role of platelets in thrombus formation drives the focus on antiplatelet therapy. The standard approach, as mentioned above, is aspirin therapy, which is simple and well-tested. However, any antiplatelet therapy balances the benefits of clot prevention with a risk of bleeding. The benefits of aspirin as a prophylactic approach against a secondary event (*e.g.*, against

#### **Arachidonic Acid**



**Figure 2.** COX-1 and COX-2 convert arachidonic acid to the intermediate  $PGH_2$ , which is then metabolized by specific distal enzymes to produce the different PGs. For example, TXA $_2$  synthase hydrolyzes  $PGH_2$  to give TXA $_2$ . TXA $_2$  is rapidly hydrolyzed non-enzymatically to TXB $_2$ .

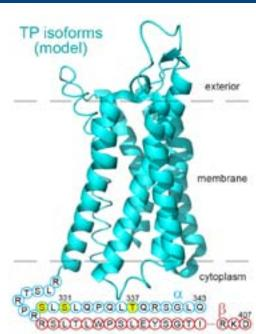


Figure 3. A model of the TP receptor isoforms, showing the critical region that differs between TP $\alpha$  (cyan) and TP $\beta$  (red). Key residues that are phosphorylated on TP $\alpha$  are highlighted in yellow.

recurrence of infarct) clearly outweighs the bleeding risk. Unfortunately, in trials of primary prevention, the benefits of this COX inhibitor only marginally outweigh the risks. Similar problems exist for flavonoids (*e.g.*, apigenin, luteolin, quercetin), which have antiplatelet effects through inhibition of the COX pathway.

TX blockers, blocking either  $TXA_2$  synthase activity or TP receptors, allow the biosynthesis of other COX pathway products, most notably  $PGI_2$ . Moreover,  $TXA_2$  synthase inhibitors prevent the generation of  $TXA_2$  in sources other than platelets. Potent inhibitors of  $TXA_2$  synthase, TP receptors, as well as dual inhibitors, are available from Cayman (Table 1). Similar compounds are being tested or used as antithrombotic agents in the treatment of peripheral arterial disease (PAD), PAD with type 2 diabetes, and stroke.

Inhibitors and Antagonists			
Item No.	Product Name		
70540	Furegrelate (sodium salt)	Inhibitor of human platelet TXA $_2$ synthase (IC $_{50}$ = 15 nM)	
70515	Ozagrel	Selective inhibitor of TXA <sub>2</sub> synthase ( $IC_{50} = 11 \text{ nM}$ )	
19025	SQ 29,548	Potent antagonist of the TP receptor: K <sub>i</sub> of 4.1 nM	
10011562	L-655,240	Potent antagonist of the TP receptor <i>in vitro</i> (IC <sub>50</sub> = 7 nM)	
10010396	CAY10535	TP receptor antagonist, with selectivity for TP $\beta$ (IC $_{50}$ $=$ 99 nM) over TP $\alpha$ (IC $_{50}$ $=$ 1,970 nM)	
10155	BM 567	A dual acting antithrombogenic agent, acting as an inhibitor of TXA $_2$ synthase (IC $_{50}=12$ nM) and as an antagonist of the TP receptor (IC $_{50}=1.1$ nM)	
For a full product listing, please visit www.caymanchem.com			

#### References

- 1. Davì, G., Catalano, I., Averna, M., et al. N. Engl. J. Med. 322, 1769-1774 (1990).
- 2. Coyle, A.T. and Kinsella, B.T. *Biochem. Pharmacol.* **71,** 1308-1323 (2006).
- 3. Gannon, A.M.M., Turner, E.C., Reid, H.M., et al. J. Mol. Biol. **3394**, 29-45 (2009).
- Yamamoto, S., Yan, F., Zhou, H., et al. Biochem. Pharmacol. 64, 375-383 (2002).
   Reid, H.M. and Kinsella, B.T. J. Biol. Chem. 278(51), 51190-51202 (2003).
- 6. Kelley-Hickie, L.P., O'Keeffe, M.B., Reid, H.M., et al. Biochem. Biophys. Acta 1773(6), 970-989 (2007).
- 7. Kelley-Hickie, L.P. and Kinsella, B.T. Br. J. Pharmacol. 142, 203-221 (2004).
- 8. Nakahata, N. Pharmacol. Ther. 118, 18-35 (2008).
- 9. Coccheri, S. Drugs **70(7)**, 887-908 (2010)

#### Hormone Sensitive Lipase (human recombinant)

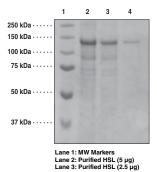
10664

HSL M<sub>r</sub>: 118.3 kDa Purity: ≥80%

Supplied in: 50 mM sodium phosphate, pH 7.2, containing 100 mM sodium chloride and 20% glycerol

Source: Recombinant N-terminal His-tagged protein, expressed in Sf21 insect cells using a baculovirus expression system

25 µg 50 µg 100 ua



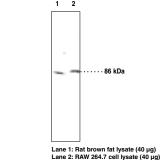
#### Hormone Sensitive Lipase Polyclonal Antibody

10006371

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: synthetic peptide from human HSL amino acids 731-741 Host: rabbit • Cross Reactivity: (+) human, murine, and rat HSL • Application(s): WB • HSL catalyzes the hydrolysis of tri-, di-, and monoacylglycerols, as well as cholesterol esters.

500 µl



#### 13(S)-HOTrE 39620

[87984-82-5]

MF:  $C_{19}H_{30}O_3$  FW: 294.4 Purity: ≥98%

A solution in ethanol **Stability:** ≥2 years at -20°C

Summary: 15-LO product of linolenic acid; detected in cell membranes and as the cholesteryl ester associated with atherosclerotic lesions

100 µg 500 µg 1 mg

5 mg

10699

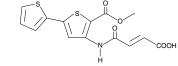
## HTS 01037

[682741-29-3]

**MF:**  $C_{14}H_{11}NO_{5}S_{2}$  **FW:** 337.4 **Purity:** ≥98% A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A high-affinity ligand of FABP4 ( $K_i = 0.67 \mu M$ ) that has been shown to antagonize the protein-protein interaction of FABP4 with hormone sensitive lipase in cultured C8PA lipocytes; it presumably competes with fatty acids for functional binding in the ligand-binding cavity of FABP4

1 mg 5 mg 10 ma 25 ma



I-BOP 19600

[128719-90-4]

**MF:**  $C_{23}H_{29}IO_5$  **FW:** 512.4 **Purity:**  $\geq$ 95%\*

A solution in ethanol **Stability:** ≥2 years at -20°C

Summary: A potent TP receptor agonist; induces human platelet aggregation with EC<sub>50</sub> values of 0.34 and 0.17 nM at pH 7.4 and pH 6.0, respectively

500 µg 1 mg 5 mg

18215

lloprost

[78919-13-8] Ciloprost

MF:  $C_{22}H_{32}O_4$  FW: 360.5 Purity:  $\geq 97\%$ 

A solution in methyl acetate **Stability:** ≥1 year at -20°C

Summary: A second generation structural analog of prostacyclin; binds with equal affinity to the human recombinant IP and EP<sub>1</sub> receptors (K<sub>i</sub> = 11 nM)

500 µg 1 mg 5 mg 10 mg

#### IP Receptor (human) Polyclonal Antibody 10005518

PGI<sub>2</sub> Receptor, PGI<sub>2</sub> Receptor, Prostacyclin Receptor Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human IP receptor N-terminal amino acids 1-16 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat IP receptor • Application(s): WB • The PGI2 receptor participates in signal transduction of the pain response, cardioprotection, and inflammation.

500 µl

Also Available:

IP Receptor (human) Blocking Peptide (10005519)

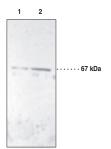
#### IP Receptor (murine) Polyclonal Antibody 160070

PGI<sub>2</sub> Receptor, PGI<sub>2</sub> Receptor, Prostacyclin Receptor

Peptide affinity-purified IgG **Stability:** ≥2 years at -20°C Summary: Antigen: murine IP receptor N-terminal amino acids 3-16 •

Host: rabbit • Cross Reactivity: (+) murine and rat IP receptor; (-) human IP receptor • Application(s): WB • The IP receptor mediates the vasorelaxation and anti-platelet aggregation effects of prostacyclin (PGI<sub>2</sub>)

500 µl



• Also Available: IP Receptor (murine) Blocking Peptide (360070)

\*All 5-cis 2-series PGs (those containing a 5,6-double bond) will contain a small amount of the 5-trans isomer. This isomer is generally undetectable using normal phase silica columns and plates, but may be resolved using RP-HPLC. The purity for all such 2-series PGs excludes the 1-3% trans isomer which will generally be present.

IP and TP Receptors Agonists and Antagonists						
Item No.	Product Name	Target	Mode of Action	Effective Concentration	Sizes	
18230	Beraprost (sodium salt)	IP	Agonist	Doses of 20-100 μg are effective in humans	1 mg • 5 mg • 10 mg • 50 mg	
10155	BM 567	TP	Antagonist	$IC_{50} = 1.1 \text{nM}$	1 mg • 5 mg • 10 mg • 50 mg	
18210	Carbaprostacyclin	IP	Agonist	$ED_{50} = 47$ nM for inhibition of platelet aggregation	500 μg • 1 mg • 5 mg • 10 mg	
19010	Carbocyclic Thromboxane A <sub>2</sub>	TP	Agonist	1 nM effectively constricts cat coronary arteries	100 μg • 500 μg • 1 mg • 5 mg	
10005186	CAY10441	IP	Antagonist	$K_i = 1.5 \text{ nM}$	1 mg • 5 mg • 10 mg • 25 mg	
10005913	CAY10449	IP	Antagonist	K <sub>i</sub> = 3 nM	1 mg • 5 mg • 10 mg • 25 mg	
10010396	CAY10535	TP	Antagonist	$IC_{50} = 99 \text{ nM (TP}\beta); IC_{50} = 1,970 \text{ nM (TP}\alpha)$	1 mg • 5 mg • 10 mg • 50 mg	
10012447	CG 4305	IP	Agonist	EC <sub>25</sub> = 25 nM for inhibition of platelet aggregation	100 μg • 500 μg • 1 mg • 5 mg	
18216	Ciprostene (calcium salt)	IP	Agonist	${\rm ID}_{\rm 50} =$ 60 ng/ml for inhibition of ADP-induced platelet aggregation <i>in vitro</i>	1 mg • 5 mg • 10 mg • 50 mg	
19600	I-BOP	TP	Agonist	EC <sub>50</sub> = 0.34 nM (pH 7.4); 0.174 nM (pH 6.0)	100 μg • 500 μg • 1 mg • 5 mg	
18215	lloprost	IP	Agonist	$K_i = 11 \text{ nM}$ (binds with equal affinity to EP <sub>1</sub> receptor)	500 μg • 1 mg • 5 mg • 10 mg	
19021	I-SAP	TP	Antagonist	$K_d = 0.5$ nM (human platelets)	100 μg • 500 μg • 1 mg • 5 mg	
10011562	L-655,240	TP	Antagonist	$IC_{50} = 7 \text{ nM}$	500 μg • 1 mg • 5 mg • 10 mg	
10010412	mRE-269	IP	Agonist	$K_i = 20 \text{ nM}$	1 mg • 5 mg • 10 mg • 50 mg	
19020	Pinane Thromboxane A <sub>2</sub>	TP	Antagonist	$IC_{50} = 2~\mu\text{M}$ for inhibition of U-46619-induced aggregation of human platelets	500 μg • 1 mg • 5 mg • 10 mg	
18220	Prostaglandin I <sub>2</sub> (sodium salt)	IP	Agonist	$IC_{50} = 5$ nM for inhibition of human platelet aggregation	1 mg • 5 mg • 10 mg • 25 mg	
19025	SQ 29,548	TP	Antagonist	$K_i = 4.1  \text{nM}$	1 mg • 5 mg • 10 mg • 100 mg	
10011348	Taprostene	IP	Agonist		500 μg • 1 mg • 5 mg • 10 mg	
16440	U-44069	TP	Agonist	$EC_{50} = 3 \mu M$ for platelet aggregation	1 mg • 5 mg • 10 mg • 50 mg	
16450	U-46619	TP	Agonist	EC <sub>50</sub> = 82 nM for human platelet aggregation	1 mg • 5 mg • 10 mg • 50 mg	

19021

[133538-58-6] Iodophenyl sulfonyl amino pinane TXA2 MF:  $C_{22}H_{30}INO_4S$  FW: 531.4 Purity:  $\geq 98\%^*$ 

A solution in ethanol **Stability:** ≥2 years at -20°C

Summary: A high-affinity TP receptor antagonist; produces platelet shape change, but not aggregation, with an EC<sub>50</sub> value of 9.7 nM

100 µg 500 µg 1 mg 5 mg

I-SAP

#### KB2115 10011054

[355129-15-6]

**MF:**  $C_{19}H_{17}Br_{2}NO_{5}$  **FW:** 487.1 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A synthetic thyroid hormone mimetic that lowers total and LDL cholesterol up to 40% when administered to humans at a dose of 50-200 µg once daily for 14 days; stimulates cholesterol catabolism to bile acids without affecting cholesterol synthesis

500 µg 1 mg 5 mg 10 mg

#### KDdiA-PC

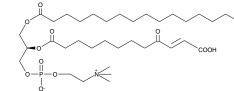
62935

[439904-34-4]

MF:  $C_{36}H_{66}NO_{11}P$  FW: 719.9 Purity:  $\geq$ 98% A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A phosphatidylcholine species containing a fragmented, oxidized shortchain fatty acid remnant at the sn-2 position; acts as a potent CD36 ligand

500 µg 1 mg 5 mg 10 mg



#### KOdiA-PC 62945

[439904-33-3]

MF:  $C_{32}H_{58}O_{11}P$  FW: 663.8 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A phosphatidylcholine species containing a fragmented, oxidized shortchain fatty acid remnant at the sn-2 position; acts as a potent CD36 ligand

500 µg 1 mg 5 mg 10 mg

#### L-655,240

10011562

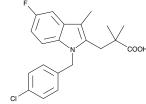
[103253-15-2]

**MF:**  $C_{21}H_{21}ClFNO_2$  **FW:** 373.9 **Purity:**  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A potent antagonist of the  $TXA_2$  receptor (IC<sub>50</sub> = 7 nM); blocks TPmediated bronchoconstriction in vivo and platelet aggregation ex vivo

500 µg 1 mg 5 mg 10 mg



#### LCAT Polyclonal Antibody

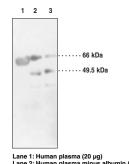
10009323

10007665

Lecithin: Cholesterol Acyltransferase, Phosphatidylcholine-Sterol O-Acyltransferase Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human LCAT amino acids 132-143 • Host: rabbit • Cross Reactivity: (+) human, murine, porcine, and bovine LCAT • Application(s): WB • LCAT catalyzes the fatty acid transfer from the sn-2 position of phosphatidylcholine (lecithin) to cholesterol and to a lesser degree to other acceptor molecules. This enzyme is critical to the process of reverse cholesterol transport or movement of cholesterol esters into HDL paricles from cells.





Also Available: LCAT Blocking Peptide (10009324)

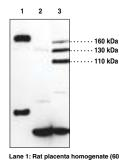
#### LDL Receptor Polyclonal Antibody

LDLR

Affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: Synthetic peptide from murine LDL receptor amino acids 499-511 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat LDL receptor Application(s): ICC and WB • The LDLRs are cell surface glycoproteins that scavenge LDL from the blood and regulate plasma LDL cholesterol.

500 ul



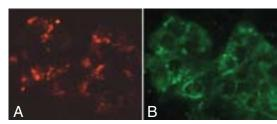
#### LDL Uptake Cell-Based Assay Kit

10011125

**Stability:** ≥6 months at 4°C

Summary: LDL uptake and its regulation are important therapeutic targets for atherosclerosis and related diseases. Cayman's LDL Uptake Cell-Based Assay employs a preparation of human LDL conjugated to Dylight<sup>TM</sup> 549 as a fluorescent probe for detection of LDL uptake into cultured cells. A LDL receptor-specific antibody and a Dylight<sup>TM</sup> 488-conjugated secondary antibody are included in the kit for identifying the distribution of LDL receptors.

1 ea



LDL Uptake in HepG2 cells. HepG2 cells were cultured for two days and then treated with 32 µM EGCG overnight. Panel A: LDL-DyLight Ti 549 taken into cells appear in red. Panel B: LDL receptors in green show a distribution pattern that matches cells in Panel A containing

#### Leukotriene A, Hydrolase (human recombinant)

 $LTA_{A}H$ M<sub>r</sub>: 69.3 kDa Purity: ≥90%

Supplied in: 100 mM Tris, pH 8.0, containing 100 mM potassium chloride and 20% glycerol Source: Recombinant C-terminal His-tagged enzyme expressed in E. coli • LTA4 hydrolase is a bifunctional zinc metalloenzyme that converts LTA4 into LTB<sub>4</sub>.

25 µg 50 µg 100 µg

• Also Available: LTA, Hydrolase Polyclonal Antibody (160250)

#### Linolein Hydroperoxides

89430

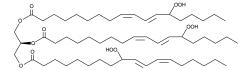
10007817

**Purity:** ≥98% (A mixture of 132 isomers)

A solution in ethanol **Stability:** ≥2 years at -80°C

Summary: A mixture of 132 possible isomers of mono-, di-, and tri-hydroperoxides produced from the autoxidation of trilinolein; may contribute to the pathophysiology of atherosclerosis when present in the circulation

500 µg 1 mg 5 mg 50 ma

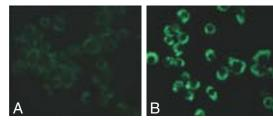


#### Lipid Droplets Fluorescence Assay Kit

**Stability:** ≥1 year at -20°C

Summary: Lipid droplets are a fundamental component of intracellular lipid homeostasis in all cell types and they provide a rapidly mobilized lipid source for many important biological processes. Cayman's Lipid Droplets Fluorescence Assay can be used to study regulators of lipid droplet biogenesis. The main advantage of this assay is that the green fluorescence of Nile Red is both very sensitive and specific for lipid droplet detection.

480 tests



Oleic Acid dramatically induces lipid droplet accumulation in neuro-2a cells. Neuro-2a cells were treated with vehicle (control, Panel A) or 400 µM oleic acid (treated, Panel B). In cells cultured with MEM medium only, there are few cells with tiny lipid droplets (Panel A). When cells were fed with 400  $\mu M$  oleic acid , most of cells accumulated numerous large lipid droplets (Panel B).

#### 5-Lipoxygenase (human recombinant)

Mr: 78 kDa Purity: Cell lysate 16,000 x g supernatant

Supplied in: 100 mM Tris, pH 8.0, containing 5 mM EGTA

**Source:** Recombinant enzyme isolated from a Baculovirus overexpression system in Sf21 cells

500 units

1 Kunit

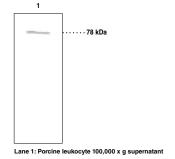
2.5 Kunit 5 Kunit

#### 5-Lipoxygenase Polyclonal Antibody

Peptide affinity-purified IgG **Stability:** ≥2 years at -20°C

Summary: Antigen: human and rat amino acids 130-149; The sequence is 95% homologous to 5-LO from mouse and hamster • Host: rabbit • Cross Reactivity: (+) human, rat, murine, hamster, and porcine 5-LO; (-) 12-LO and 15-LO • Application(s): ICC, IHC, and WB • 5-LO catalyzes the formation of 5(S)-HpETE from arachidonic acid as well as its subsequent conversion to LTA<sub>4</sub>.

500 µl



• Also Available: 5-Lipoxygenase Blocking Peptide (360402)

#### 5-Lipoxygenase (Phospho-Ser<sup>523</sup>) Polyclonal Antibody

**Stability:** ≥1 year at -20°C

Summary: Antigen: phosphopeptide corresponding to amino acid residues surrounding the phospho-Ser<sup>523</sup> of human 5-LO • Host: rabbit • Cross Reactivity: (+) human, rat, and non-human primate 5-LO • Application: WB • The 5-LO enzyme plays a key role in regulating the production of LTs. Its activity is regulated by PKA phosphorylation at serine-523.

1 ea

#### 12-Lipoxygenase (murine leukocyte) Polyclonal Antiserum

160304

Lyophilized antiserum **Stability:** ≥2 years at -20°C

Summary: Antigen: recombinant murine leukocyte 12-LO • Host: rabbit • Cross Reactivity: (+) murine, porcine, and human leukocyte 12-LO and rabbit reticulocyte 15-LO • Application(s): WB

500001

60402

160402

10007820

#### 12-Lipoxygenase (platelet-type, murine recombinant)

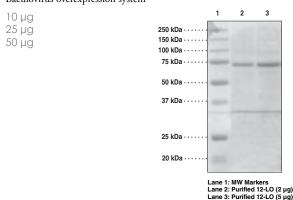
10341

12-LO

M<sub>.</sub>: 76.4 kDa Purity: ≥85%

Supplied in: 50 mM sodium phosphate, pH 7.2, containing 100 mM sodium chloride, 20 µM ferrous chloride, and 30% glycerol

**Source:** Recombinant N-terminal His-tagged protein, expressed in Sf21 cells using a Baculovirus overexpression system



#### 15-Lipoxyaenase-1 (rabbit) Polyclonal Antiserum

160704

16833

Lyophilized antiserum **Stability:** ≥3 years at -20°C

Summary: Antigen: rabbit reticulocyte 15-LO • Host: sheep • Cross Reactivity: (+) rabbit, human, and murine 15-LO-1 and 12-LO (porcine leukocyte) • Application(s): WB • 15-LO-1 catalyzes the formation of 15(S)-HETE and 13(S)-HODE from arachidonic acid and linoleic acid, respectively.

500 ul

#### Lisinopril

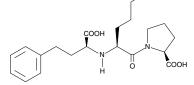
[76547-98-3] MK 521, MK 522

and canine, respectively

MF:  $C_{21}H_{21}N_{2}O_{5}$  FW: 405.5 Purity:  $\geq$ 98% A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A potent ACE inhibitor that blocks the formation of angiotensin I with an IC<sub>50</sub> value of 1.2 nM (ovine) and displays ID<sub>50</sub> values of 2.3 and 6.5 μg/kg in rat

100 mg 500 mg 1 g 5 g



<sup>•</sup> Also Available: LDL Receptor Blocking Peptide (10007672)

26 | Cayman Chemical | caymanchem.com Cayman Chemical

# Think of Cayman as Epigenetics Central

In 1998, Vincent Felitti and colleagues asked over 13,000 adults to complete a questionnaire addressing their personal adverse childhood experiences (ACE), experiences that included psychological, physical, or sexual abuses as well as indicators of family dysfunction. The questionnaire responses were then compared to measures of each individual's adult risk behavior, health status, and disease. Remarkably, this first "ACE" study found a "strong graded relationship between the breadth of exposure to abuse or household dysfunction during childhood and multiple risk factors for several of the leading causes of death in adults." That is, the greater the ACE score, the more likely the individual would, as an adult, smoke or use illicit drugs, suffer from alcoholism or depression, or attempt suicide. Moreover, increasing ACE scores correlated with increased prevalence and risk for ischemic heart disease, cancer, stroke, and diabetes. This landmark study has since been corroborated by both retrospective and prospective studies that link adverse psychosocial experiences during childhood with an increased risk for age-related diseases much later in life.<sup>2,3</sup>

In 1999, researchers looking for genes that were linked to cancer began to discover gene products that acted as methyltransferases or demethylases, acetyltransferases or deacetylases. In the intervening years, an explosion of research has developed our understanding of epigenetics, the persistent modification of nucleosomes or genetic material to alter gene expression without changing DNA sequence or copy number. For example, the methylation of histones, which commonly alters gene expression, can persist through mitosis and contribute to cell differentiation, organogenesis, and cancer (see related story, page 12).

Today, the epidemiological ACE studies converge with epigenetic models to suggest new explanations for psychosocial, as well as physiological, abnormalities in adults. For example, childhood experiences might produce critical epigenetic changes that persist into adulthood and contribute to depression. Experimentally, differences in maternal care of pups in rats alter DNA methylation and histone acetylation, affecting neural development.<sup>4</sup> Such experiments suggest new physiological mechanisms for certain diseases, like adult depression, which in turn offers hope for new therapeutic interventions. Cayman anticipates a growing interest in epigenetic research and the search for new compounds that modulate epigenetic pathways. As a result, we offer multiple avenues for epigenetic discovery.

#### **Histone Acetylation**

The term 'epigenetics' refers, in its broadest sense, to persistent changes in gene expression, exclusive of changes in DNA sequence, DNA copy number, or changes due directly to transcription factors. Much contemporary excitement focuses on the nucleosome. An old model presented this structure as a spool of histones with a thread of DNA wrapped around it, serving a physical mechanism for organizing the extensive genetic material within the nucleus (Figure 1). Now, it is clear that the nucleosome is a sophisticated device for tying down or feeding out genetic material and regulating transcription. One way to understand this 'device' involves considering the positive and negative charges distributed along DNA and histone components. From this viewpoint, the DNA backbone is a string of negatively charged phosphate groups. Histones are rich in basic residues, particularly at the amino termini, which contain regularly-spaced lysine residues (see related article, page 12). These positive charges set up an electrostatic potential between the DNA and histones, defining a physical interaction that binds the thread to the spool. That is, unmodified lysines can clamp down on DNA, preventing its access to promoters. The positive charges can be neutralized by methylation or reduced by acetylation. In addition, serines, threonines, and tyrosines can be phosphorylated, adding negative charges to specific sites on the histones. In Figure 1, lysine residues are colored red, while serines known to be phosphorylated are black.

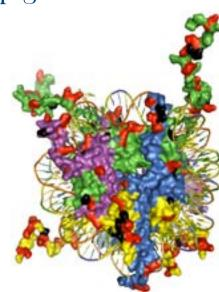


Figure 1. Nucleosomal structure, displaying the histones H2A (blue), H2B (yellow), H3 (green), and H4 (purple) in surface mode. Lysines known to be modified are colored red;

Interestingly, the modification of histones can range from crude, with multiple sites being altered en masse, to elegant and complex, with changes in individual sites controlled by sophisticated protein complexes under

A large number of acetyltransferases add an acetyl group (CH3CO) to proteins, including histones. The acetylation of histones typically relaxes chromatin conformation, facilitates the recruitment of effector proteins, and increases transcription. These histone acetyltransferases (HAT) are also known as 'lysine (K) acetyltransferases' (KAT), since they can target proteins other than histones. Acetyl groups are removed by histone deacetylases (HDAC, a.k.a. KDAC). The human class I HDACs are homologous to the yeast enzyme Rpd3 and include HDAC1, 2, 3, and 8. Class II HDACs are homologous to yeast HDA1 and are divided into class IIa (HDAC4, 5, 7, 9) and class IIb (HDAC6, 10) based on structure. The human class III HDACs include the sirtuin (SIRT) family of NAD+-dependent protein deacetylases. The novel HDAC11 has a distinct structure and is a class IV HDAC. The HDACs often participate in the formation of transcriptional complexes, inducing chromatin compaction through histone deacetylation and silencing gene expression.

HDACs, including SIRTs, have diverse roles in development and pathogenesis. 5-7 Most prominently, HDAC inhibitors affect inflammatory signaling. In humans, they inhibit NF-κB activation, while in mice and rats, they reduce the expression of tumor necrosis factor- $\alpha$  (TNF- $\alpha$ ), a potent activator of NF-kB. HDAC inhibitors also suppress pro-inflammatory cytokine expression and block T-cell infiltration. HDACs are also necessary for normal cardiovascular development and endothelial functioning, with class I HDACs playing key roles in heart and cardiomyocyte development and HDAC7 contributing to endothelium integrity. SIRTs have principle roles in metabolic and inflammatory pathways, with SIRT1 and SIRT7 being important for cardiomyocyte survival.8,9

Cayman Chemical is a leader in the development of research tools for epigenetics research. The Cayman Protein Core produces pure, functional recombinant proteins and protein complexes. Many recombinant HDACs and SIRTs, as well as several histones, are currently available (Table 1). In addition, the Cayman Assay Development Team creates versatile and dependable assay kits. We also offer antibodies as well as a broad selection of pure and potent research chemicals.

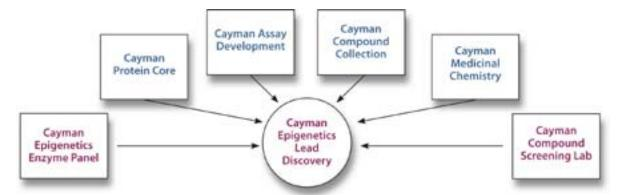


Figure 2. Cayman has historical strengths in recombinant protein production, assay development, compound synthesis, and medicinal chemistry. These are combined with a growing Epigenetics Enzyme Panel and a high-throughput Compound Screening Lab to drive Epigenetics Lead Discovery in our new facility.

Cardiovascular complications are the major cause of morbidity and mortality in the diabetic population. Exposure to high glucose is a major factor leading to these complications. Treatment is confounded by a well-documented phenomenon called "metabolic memory", which refers to the ability of certain organs to 'remember' their early glycemic environment and 'behave' accordingly.<sup>10</sup> For example, diabetic dogs switched from poor glucose control (no insulin treatment) to good control at two months, when examined at five years after the start of the study, suffered little retinopathy. Diabetic animals switched to good control at 2.5 years had a similar incidence of retinopathy as their control counterparts who received poor glucose control throughout the 5 year study. Many studies like this, both in the lab and the clinic, reflect the challenges of preventing cardiovascular complications through glucose control in diabetes.

Recent studies suggest that methyltransferases and demethylases contribute to inflammation during hyperglycemia that intertwines with subsequent vascular damage during diabetes. 11-13 In this model, acute hyperglycemia increases monomethylation of lysine 4 on histone 3 (H3K4) and suppresses methylation of H3K9 by altering the activities of the methyltransferases SET7/9 and SUV39H1, respectively. These changes lead to persistent activation of NF-kB through the enhanced recruitment of p65, upregulating pro-inflammatory pathways that drive cardiovascular complications. Most importantly, for these changes to lie at the heart of metabolic memory, they must be made at an early, impressionable stage and persist through multiple cell divisions over many years. It remains to be clarified how the methylation status of H3K4 and H3K9 can be faithfully preserved through mitosis and over time. Moreover, these changes must be consistently and specifically tied to NF-κB expression.

SET7/9 is also known as KMT7, for lysine methyltransferase 7. Similarly, LSD1 is also known as KDM1, for lysine demethylase 1. These newer names reflect a growing appreciation for the fact that both methyltransferases and demethylases target many proteins in addition to histones. For example, methylation of p53 by SET7/9 increases p53 activity. In fact, SET7/9 has numerous non-histone protein targets. 14 Is it possible that acute hyperglycemia at a formative age produces an epigenetic change on the expression of SET7/9 itself, perhaps through DNA methylation, and that metabolic memory involves SET7/9-dependent methylation of multiple proteins, including histones linked to NF-κB?

These and other questions involving methyltransferases and demethylases can be addressed using reagents available from Cayman. Cayman has created a full-service Epigenetics Lead Discovery facility that will identify small-molecule modulators of epigenetic target enzymes (Figure 2). This facility houses Cayman's new Compound Screening Lab, which will enable high-throughput screening of our in-house compound library (a 14,400 compound Maybridge HitFinder collection) or user-supplied collections against an individual target or a panel of epigenetic enzymes. To support these endeavors, additional recombinant proteins and assays are in development, with new products being added even as this article goes to press. Visit caymanchem.com to discover what's new today!

Histone acetylation/deacetylation proteins, kits, antibodies, and reagents			
Category	Product (Item No.)		
HDACs (human recombinant)	HDAC1 (1009231), HDAC2 (10009377), HDAC3/NCOR2 (10009232), HDAC4 (10009652), HDAC5 (10009379), HDAC6 (10009465), HDAC8 (19380), HDAC9 (10009466)		
SIRTs (human recombinant)	SIRT1 (10011190), SIRT2 (10011191), SIRT3 (10011194), SIRT4 (10317), SIRT5 (10318), SIRT6 (10315), SIRT7 (10316)		
Histones and Acetyltransferases (recombinant)	Histone H2B (Xenopus) (10262), Histone H3 (G103A substitution) (10263), Histone H4 (10264), pCAF Histone Acetyltransferase (10009115)		
Assay Kits	HAT Inhibitor Screening Assay Kit (10006515), HDAC Activity Assay Kit (10011563), HDAC Cell-Based Activity Assay Kit (600150), HDAC1 Inhibitor Screening Assay Kit (10011564), HDAC8 Inhibitor Screening Assay Kit (700230), SIRT1 Direct Fluorescence Screening Assay Kit (10010401), SIRT1 FRET-Based Screening Assay Kit (10010991), SIRT2 Direct Fluorescence Screening Assay Kit (700280), SIRT3 Direct Fluorescence Screening Assay Kit (7001566), SIRT6 Direct Fluorescence Screening Assay Kit (700290)		
Antibodies	HDAC1 (13491), HDAC3 (13493), HDAC4 (13494), HDAC6 (13499), HDAC7 (Phospho-Ser <sup>155</sup> ) (13500), HDAC11 (13504), TIP60 (13789), Acetyl Lysine (Clone 7F8)(10010567), SIRT7 (13477)		
Key Reagents	CAY 10398 (89740), CAY 10591 (10009797), CAY10603 (13146), CBHA (13172), Chidamide (13686), coumarin-SAHA (10671), EX-527 (10009798), (S)-HDAC-42 (13277), HNHA (13295), M 344 (13174), 4-iodo-SAHA (10495), MS-275 (13284), trans-Resveratrol (70675), SAHA (10009929), Salermide (13178), Scriptaid (10572), Sodium Butyrate (13121), SRT 1720 (10011020), Tenovin-1 (13085), Tenovin-6 (13086), Trichostatin A (89730), Tubastatin A (trifluoroacetate salt) (10559), UNCO224 (13631), UNCO321 (10582)		

- Felitti, V.J., Anda, R.F., Nordenberg, D., et al. Am. J. Prev. Med. 14, 245-258 (1998).
- Dong, M., Giles, W.H., Felitti, V.J., et al. Circ. J. 110, 1761-1766 (2004).
- Danese, A., Moffitt, T.E., Harrington, H., et al. Arch. Pediatr. Adolesc. Med. 163(12), 1135-1143 (2009).
- Zhang, T.-Y., Hellstrom, I.C., Bagot, R.C., et al. J. Neurosci. 30(39), 13130-13137 (2010)
- Fernandez, A.Z., Siebel, A.L., and Fl-Osta, A. Int. J. Vasc. Med. [Unpublished] (2011)
- Haberland, M., Montgomery, R.L., and Olson, E.N. Nat. Rev. Genet. 10, 32-42 (2011)
- Blum, C.A., Ellis, J.L., Loh, C., et al. J. Med. Chem. 54, 417-432 (2011).
- Sundaresan, N.R., Pillai, V.B., and Gupta, M.P. J. Mol. Cell. Cardiol. [Unpublished] (2011). Vakhrusheva, O., Smolka, C., Gajawada, P., et al. Circ. Res. 102, 703-710 (2008).
- Ceriello, A., Ihnat, M.A., and Thorpe, J.E. J. Clin. Endocrinol. Metab. 94(2), 410-415 (2009).
- I. Li, Y., Reddy, M.A., Mlao, F., et al. J. Biol. Chem. 283(39), 26771-26781 (2008).
- Brasacchio, D. Okabe, I. Tikellis, C., et al. Diabetes 58, 1229-1236 (2009). 13. Siebel, A.L., Fernandez, A.Z., and El-Osta, A. *Biochem. Pharmacol.* **80,** 1853-1859 (2010).
- 13. Dhayalan, A., Kudithipudi, S., Rathert, P., et al. Chem. Biol. 18, 111-120 (2011)

#### Losartan (potassium salt)

[124750-99-8] DuP 753, MK 954

**MF:** C<sub>22</sub>H<sub>23</sub>ClN<sub>6</sub>O • K **FW:** 461.0 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An  $AT_1$  receptor antagonist ( $K_i = 5-20$  nM); attenuates vein graft atherosclerosis in rabbits and reduces arterial blood pressure in rats; controls hypertension while protecting renal function in humans

caymanchem.com

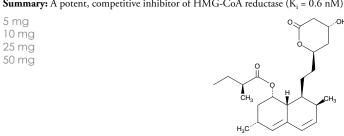
10 mg 50 mg 100 mg 500 mg

#### Lovastatin [75330-75-5]

MF:  $C_{24}H_{36}O_5$  FW: 404.5 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent, competitive inhibitor of HMG-CoA reductase  $(K_i = 0.6 \text{ nM})$ 



• Also Available: Lovastatin Hydroxy Acid (sodium salt) (10010339)

#### Luminex® Thromboxane B<sub>2</sub> Kit

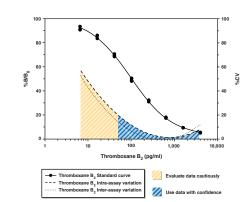
10007502

**Stability:** ≥1 year at -20°C

**Sensitivity:** 80% B/B<sub>0</sub>: ~24 pg/ml • 50% B/B<sub>0</sub>: ~100 pg/ml

Summary: Cayman's Luminex<sup>□</sup> TXB<sub>2</sub> can be used for quantification of TXB<sub>2</sub> in plasma, culture media, and other sample matrices.

1 ea



#### Luminex® 11-dehydro Thromboxane B<sub>2</sub> Kit 10010971

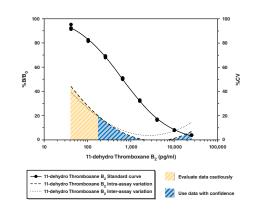
**Stability:** ≥1 year at -20°C

**Sensitivity:** 50% B/B<sub>0</sub>=650 pg/ml • 80% B/B<sub>0</sub>=130 pg/ml

Summary: 11-dehydro TXB<sub>2</sub> is a metabolite of TXB<sub>2</sub> which can be measured in urine to reflect systemic production of TXB2 in vivo. Cayman's Luminex 11-dehydro TXB2 is the first of its kind for the measurement of 11-dehydro TXB2 using Luminex<sup>□</sup> xMAP<sup>□</sup> technology.

1 ea

10006594



#### MEDICA 16

90290

#### [87272-20-6]

MF:  $C_{20}H_{38}O_4$  FW: 342.5 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥1 year at -20°C

**Summary:** A β,β'-dimethyl hexadecanedioic acid that exhibits hypolipidemic and antidiabetogenic effects

I mg	
5 mg	
10 mg	
50 mg	

# Mevastatin

10010340

10010412

[73573-88-3] Compactin, CS 500, L 637312, ML 236B, NSC 281245, Statin I MF:  $C_{23}H_{34}O_5$  FW: 390.5 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A reversible, competitive HMG-CoA reductase inhibitor  $(K_i = 1 \text{ nM})$ 

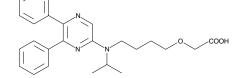


#### MRE-269 [475085-57-5]

MF:  $C_{25}H_{29}N_3O_3$  FW: 419.5 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A stable analog of prostacyclin that potently and selectively activates the IP receptor (K<sub>i</sub> = 20 nM); active form of the prodrug NS-304 that is stable in vivo, as plasma concentrations of MRE-269 remain near peak levels for more than eight hours; exhibits lower affinity for the vasoconstricting EP3 receptor than other IP receptor agonists

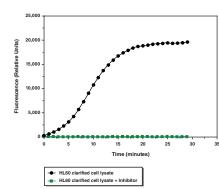


#### Myeloperoxidase Chlorination Assay Kit 10006438

**Stability:** ≥6 months at -20°C

Summary: Cayman's Myeloperoxidase Chlorination Assay provides a convenient fluorescence-based method for detecting the MPO chlorination activity in both crude cell lysates and purified enzyme preparations. The assay utilizes the non-fluorescent probe, APF, which is selectively cleaved by hypochlorite to yield the highly fluorescent compound fluorescein. The kit includes an MPO-specific inhibitor for distinguishing MPO activity from MPO-independent fluroescence.

2 x 96 wells



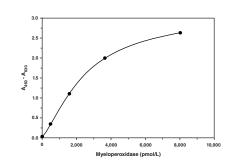
#### Myeloperoxidase (human) EIA Kit

**Stability:** ≥6 months at 4°C

Limit of Detection: 14 pg/ml

Summary: Cayman's MPO (human) EIA is an immunometric (i.e. sandwich) assay which can be used to measure MPO in plasma without prior sample purification. This assay has been tested using plasma from healthy volunteers and the results were shown to be consistent with published data.

96 wells

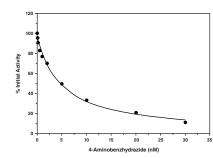


#### Myeloperoxidase Inhibitor Screenina Assav Kit

**Stability:** ≥6 months at -20°C

Summary: Cayman's MPO Inhibitor Screening Assay provides fluorescence-based methods for screening inhibitors to both the chlorination and peroxidation activities of MPO. Sufficient reagents are provided for a full 96-well plate assay of each type of activity.

2 x 96 wells

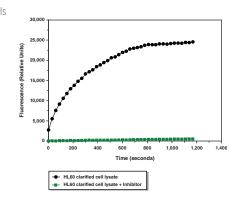


#### Myeloperoxidase Peroxidation Assay Kit

Stability: ≥6 months at -20°C

Summary: Cayman's MPO Peroxidation Assay provides a fluorescence-based method for detecting MPO peroxidase activity in both crude cell lysates and purified enzyme preparations. The MPO-catalyzed reaction between hydrogen peroxide and ADHP produces the highly fluorescent compound resorufin. The kit includes an MPO-specific inhibitor for distinguishing MPO activity from MPO-independent fluorescence.

2 x 96 wells



#### Myricetin

585001

700170

10012600

[529-44-2] Cannabiscetin, NSC 407290

MF:  $C_{15}H_{10}O_8$  FW: 318.2 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A flavonoid compound that acts as a powerful antioxidant; inhibits TBARS formation with an IC<sub>50</sub> value of 6.34 µM; blocks oxLDL uptake by U937derived macrophages at 20 µM



#### S-NEPC 10008609

MF:  $C_{16}H_{13}NO_6$  FW: 315.3 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A colorimetric substrate for measurement of sEH activity

1 mg 5 mg 10 mg 25 mg

#### eNOS (bovine recombinant)

ecNOS, NOS III M.: 135 kDa Purity: Cell lysate 100,000 x g supernatant

Supplied in: 50 mM HEPES, pH 7.4, containing 10% glycerol, 5 mM CHAPS,

Source: Recombinant enzyme isolated from a Baculovirus overexpression system in Sf9 cells

10 units

#### eNOS Polyclonal Antiserum

160880

60880

ecNOS, NOS III

Lyophilized antiserum **Stability:** ≥2 years at -20°C

Summary: Antigen: human eNOS amino acids 1186-1203 • Host: rabbit • Cross Reactivity: (+) bovine and human eNOS; (-) iNOS and nNOS • Application(s): IP and WB • eNOS catalyzes the formation of NO from L-arginine in many cells and tissues including vascular endothelium, bronchiolar epithelium, cardiac myocytes, spleen, and kidney.

500 µl

• Also Available: eNOS Blocking Peptide (360881) eNOS Electrophoresis Standard (360880)

#### NS-304 10010411

[475086-01-2] ACT-293987, Selexipag

MF:  $C_{26}H_{32}N_4O_4S$  FW: 496.6 Purity:  $\geq$ 98%

A solution in methyl acetate **Stability:** ≥1 year at -20°C

**Summary:** A long-acting and potent IP receptor agonist prodrug ( $K_i = 20 \text{ nM}$ )

5 mg 10 mg 50 mg

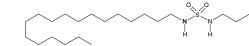
#### N-Octadecyl-N'-propyl-sulfamide 10009661

MF:  $C_{21}H_{46}N_2O_2S$  FW: 390.7 Purity:  $\geq$ 95%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A selective and potent activator of PPARa with an EC50 value of 100 nM; induces satiety, decreases food-intake, reduces body weight, and lowers plasma triglyceride concentration in free-feeding Wistar and obese Zucker (fa/fa) rats

10 mg 25 ma 50 mg



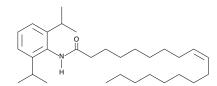
#### Oleic Acid-2,6-diisopropylanilide

[140112-65-8]

MF: C<sub>30</sub>H<sub>51</sub>NO FW: 441.7 Purity: ≥98%

A solution in methyl acetate **Stability:** ≥1 year at -20°C Summary: An inhibitor of ACAT with an IC<sub>50</sub> value of 7 nM

5 mg 10 mg 50 mg 100 ma



#### Oleoyl Ethanolamide

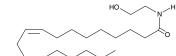
[111-58-0] OEA, Oleic Acid Ethanolamide

MF:  $C_{20}H_{30}NO_2$  FW: 325.5 Purity:  $\geq$ 98%

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: An analog of AEA found in brain tissue and chocolate; accumulates rapidly in infarcted tissue; an endogenous, potent PPAR $\alpha$  agonist (EC<sub>50</sub> = 120 nM); suppresses food intake and reduces weight gain in rats

5 ma 10 mg 50 mg 100 mg



#### Olevl Anilide

10006529

10006782

90265

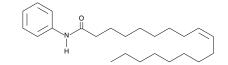
[5429-85-6] OA, Oleic Acid Anilide

**MF:** C<sub>24</sub>H<sub>39</sub>NO **FW:** 357.6 **Purity:** ≥95%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A weak ACAT inhibitor with an IC<sub>50</sub> value of 26 µM

10 ma 50 mg 100 mg



#### Oxidized Lipid HPLC Mixture

34004

Purity: ≥98% for each compound A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: Contains the free acid forms of racemic 15-HETE, 9-HODE, and 13-HODE, as well as racemic 9-HODE and 13-HODE cholesteryl esters (5 µg each)

#### oxLDL-β<sub>2</sub>GPI (human) ELISA Kit

10007893

Oxidized low-density lipoprotein- $\beta_2$  Glycoprotein I (human)

**Stability:** ≥6 months at 4°C

Summary: oxLDL is the principal form of cholesterol that accumulates in atherosclerotic lesions or plaques. Unlike native LDL, oxLDL binds to β<sub>2</sub>GPI to form oxLDL-β<sub>2</sub>GPI complexes. Stable oxLDL-β<sub>2</sub>GPI complexes are regarded as pathogenic and appear to be highly clinically relevant. Cayman's oxLDL-\(\beta\_2\)GPI (human) ELISA is an immunometric (i.e., sandwich) assay that detects the circulating oxLDL-β<sub>2</sub>GPI complex in human serum or plasma.

#### Ozagrel

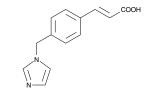
70515

[82571-53-7] OKY-046

MF:  $C_{13}H_{12}N_2O_2$  FW: 228.3 Purity:  $\geq 98\%$ A crystalline solid **Stability:** ≥1 year at -20°C

**Summary:** A selective inhibitor of TX synthase ( $IC_{50} = 11 \text{ nM}$ )

10 mg 50 mg 100 mg



#### PAF Acetylhydrolase Assay Kit

760901

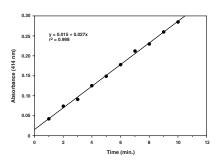
10279

Lipoprotein PLA, Lp-PLA, PAF-AH

**Stability:** ≥1 year at -20°C

Summary: PAF-AH catalyzes the hydrolysis of the potent biologically-active phospholipid PAF, generating inactive lyso-PAF. Cayman's PAF-AH Assay provides an accurate and convenient method for measurement of PAF-AH activity. The assay uses 2-thio PAF which serves as a substrate for PAF-AH. Upon hydrolysis of the acetyl thioester bond at the sn-2 position by PAF-AH, free thiols are detected using DTNB (Ellman's reagent).

96 wells



#### PAF Acetylhydrolase (human recombinant)

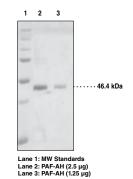
Lipoprotein PLA, Lp-PLA, PAF-AH

**M**.: 46.4 kDa **Purity:** ≥95%

Supplied in: 100 mM sodium phosphate, pH 7.2, containing, 100 mM sodium chloride and 20% glycerol

Source: Recombinant N-terminal His-tagged protein purified from E. coli

25 µg 50 µg 100 µg



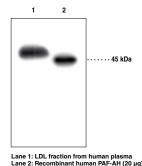
#### PAF Acetylhydrolase (human) Polyclonal Antibody

Lipoprotein PLA, Lp-PLA, PAF-AH

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human PAF-AH C-terminal amino acids 420-441; the peptide sequence used as an antigen is 64% homologous to the corresponding bovine sequence • Host: rabbit • Cross Reactivity: (+) human plasma PAF-AH; (-) murine, guinea pig, canine, and chicken PAF-AH • Application(s): WB • PAF-AH converts PAF to the biologically inactive lyso-PAF. Plasma PAF-AH is highly selective for phospholipids with very short acyl groups at the sn-2 position and is associated with lipoproteins. It has been linked to atherosclerosis and may be a positive risk factor for coronary heart disease in humans.

500 µl



• Also Available: PAF Acetylhydrolase (human) Blocking Peptide (360603) PAF Acetylhydrolase (human) Western Ready Control (10010081)

#### PAF Acetylhydrolase Inhibitor Screening Assay Kit

10004380

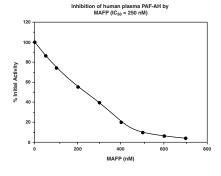
160603

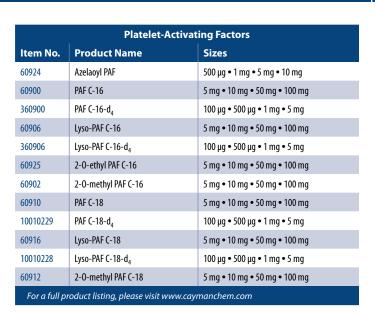
Lt-PLA, PAF-AH

**Stability:** ≥6 months at -20°C

Summary: Cayman's PAF-AH Inhibitor Screening Assay uses 2-thio PAF as a substrate for PAF-AH. Upon hydrolysis of the acetyl thioester bond at the sn-2 position by PAF-AH, free thiols are detected using DTNB (Ellman's reagent).

96 wells





#### PAF Receptor (human) Monoclonal Antibody (11A4, Clone 21)

160600

Protein-A purified IgG<sub>2a</sub> **Stability:** ≥18 months at 4°C

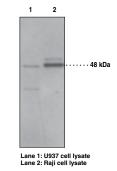
Summary: Antigen: human PAF receptor amino acids 260-269 • Host: mouse, 11A4 clone 21 • Cross Reactivity: (+) human, bovine, and porcine PAF receptor • Application(s): FC, ICC; does not work for WB • PAF is a biologically active phospholipid whose biological effects include activation of platelets, polymorphonuclear leukocytes, monocytes, and macrophages. PAF increases vascular permeability, decreases cardiac output, induces hypotension, and stimulates uterine contraction.

• Also Available: PAF Receptor Blocking Peptide (Monoclonal) (360600)

#### PAF Receptor (human) Polyclonal Antibody 160602

Peptide affinity-purified IgG **Stability:** ≥1 year at 4°C

Summary: Antigen: human PAF receptor amino acids 1-17 • Host: rabbit • Cross Reactivity: (+) human, murine, rat, and porcine PAF receptor • Application(s): FC, ICC, and WB • PAF is a biologically active phospholipid whose biological effects include activation of platelets, polymorphonuclear leukocytes, monocytes, and macrophages. PAF increases vascular permeability, decreases cardiac output, induces hypotension, and stimulates uterine contraction.



• Also Available: PAF Receptor Blocking Peptide (Polyclonal) (160604)

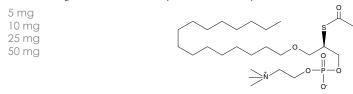
#### 2-thio-PAF 60945

[96801-55-7]

MF: C<sub>26</sub>H<sub>54</sub>NO<sub>6</sub>PS FW: 539.8 Purity: ≥98%

A solution in ethanol **Stability:** ≥1 year at -80°C

Summary: A PAF receptor agonist with potency comparable to PAF C-18 and PAF C-16; chromogenic substrate used in Cayman's PAF-AH Assay Kit (Item No. 760901)



#### Palmitoleic Acid

10009871

[373-49-9] 9-cis-Hexadecenoic Acid, Palmitoleate, n-7 Palmitoleate, cis-Palmitoleic Acid **MF:**  $C_{16}H_{30}O_{2}$  **FW:** 254.4 **Purity:** ≥99%

A neat oil **Stability:** ≥1 year at -20°C

Summary: An ω-7 monounsaturated fatty acid that is a common constituent of the triglycerides of human adipose tissue; raises LDL cholesterol and lowers HDL cholesterol much like that of a saturated fatty acid, even when dietary intake of cholesterol is maintained at a low level



#### Pargyline (hydrochloride)

10007852

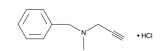
[306-07-0]

MF: C<sub>11</sub>H<sub>13</sub>N • HCl FW: 195.7 Purity: ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An irreversible inhibitor of monoamine oxidase (MAO) that is used clinically to treat moderate hypertension

100 mg 500 mg 1 g 5 g



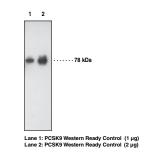
#### PCSK9 Monoclonal Antibody (Clone 15A6)

NARC-1, Proprotein Convertase Subtilisin Kexin 9

Lyophilized IgG<sub>1</sub> **Stability:** ≥1 year at -20°C

Summary: Antigen: purified human recombinant PCSK9 • Host: mouse, clone 15A6 • Isotype: IgG<sub>1</sub> • Cross Reactivity: (+) human recombinant PCSK9 • Application(s): PCSK9 ligand blotting (nondenaturing conditions) and WB • PCSK9 is a member of the subtilisin serine protease family with an important role in lipoprotein metabolism. Gain-of-function mutations in the PCSK9 gene are associated with autosomal dominant hypercholesterolemia which is characterized by an increase in LDL cholesterol levels.

1 ea



#### PCSK9 Polyclonal Antibody

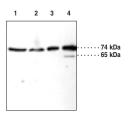
10240

NARC-1, Proprotein Convertase Subtilisin Kexin 9

Protein A-purified IgG Stability: ≥1 year at -20°C

Summary: Antigen: human recombinant PCSK9 • Host: rabbit · Cross Reactivity: (+) human, rat, and murine PCSK9 • Application(s): ICC and WB • PCSK9 is a member of the subtilisin serine protease family with an important role in lipoprotein metabolism. Gain-of-function mutations in the PCSK9 gene are associated with autosomal dominant hypercholesterolemia which is characterized by an increase in LDL cholesterol levels.

500 µl



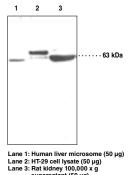
Also Available: PCSK9 Western Ready Control (10009567)

#### PCSK9 (human) Polyclonal Antibody 10007185

NARC-1, Proprotein Convertase Subtilisin Kexin 9

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: synthetic peptide from human PCSK9 amino acids 490-502 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat PCSK9 • Application(s): ICC and WB • PCSK9 is a member of the subtilisin serine protease family with an important role in lipoprotein metabolism. Several gain-of-function mutations in the PCSK9 gene are associated with hypercholesterolemia which is characterized by an increase in LDL cholesterol levels.



• Also Available: PCSK9 (human) Blocking Peptide (10007186)

#### 10008811 PCSK9 (murine) Polyclonal Antibody

NARC-1, Proprotein Convertase Subtilisin Kexin 9

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: synthetic peptide from mouse PCSK9 amino acids 152-163 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat PCSK9 • Applications: ICC and WB • PCSK9 is a member of the subtilisin serine protease family with an important role in lipoprotein metabolism. Several gain of function mutations in the PCSK9 gene are associated with hypercholesterolemia which is characterized by an increase in LDL cholesterol levels.

500 ul

• Also Available: PCSK9 (murine) Blocking Peptide (10009581)

#### **PGPC** 10044

[89947-79-5]

**MF:** C<sub>29</sub>H<sub>56</sub>NO<sub>10</sub>P **FW:** 609.7 **Purity:** ≥98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A predominant low molecular weight species of oxidized LDL; induces the expression of both E-selectin and VCAM-1, and increases endothelial cell binding by both neutrophils and monocytes

#### 10009134 **PHOME**

NOTE: This substrate should only be used with the pure EH

MF:  $C_{23}H_{19}NO_4$  FW: 373.4 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A fluorogenic, sensitive substrate for human sEH which displays good aqueous stability and solubility making it ideal for high throughput screening programs

#### Phosphatidylcholine Assay Kit

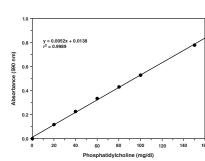
1 mg

5 mg

**Stability:** ≥6 months at -20°C

Summary: Cayman's PC Assay provides a specific, sensitive, and convenient method for quantifying PC in plasma or serum. In this assay, PC-specific PLD hydrolyzes PC to choline and phosphatidic acid. A coupled enzymatic reaction system generates a blue dye with maximum absorption at 595 nm.

96 wells



#### Pinane Thromboxane A<sub>2</sub>

[71111-01-8] PTA<sub>2</sub>

MF:  $C_{24}H_{40}O_3$  FW: 376.6 Purity:  $\geq 98\%^*$ 

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: A stable analog of TXA2 that acts as a TP receptor antagonist and inhibitor of TX synthase

#### POV-PC

[121324-31-0] 2-(5-oxovaleryl) Phosphatidylcholine

MF:  $C_{29}H_{56}NO_9P$  FW: 593.7 Purity:  $\geq$ 98%

A solution in ethanol **Stability:** ≥1 year at -20°C

Summary: One of the oxLDL species derived from 2-arachidonoyl or eicosapentanoyl phospholipids; confers CD36 scavenger receptor binding affinity of oxLDL

#### PPARα LBD (human recombinant) 10009088

PPARα Ligand Binding Domain

M\_: -34 kDa Purity: ≥90%

Supplied in: 50 mM sodium phosphate, pH 7.2, containing 20% glycerol, 100 mM sodium chloride, and 1 mM DTT

Source: Recombinant His-tagged protein purified from E. coli

50 µg 100 µg

### PPARα Polyclonal Antibody

101710

10007451

Peroxisome Proliferator-activated Receptor α

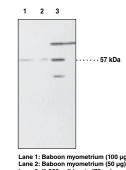
Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human, mouse, and rat amino acids 22-36 • Host: rabbit • Cross Reactivity: (+) human, murine, rat, ovine, and porcine PPARα; (-) PPARγ • Application(s): WB • PPARα is a ligand-activated transcription factor involved in the regulation of lipid homeostasis.

500 µl

10009926

19020



• Also Available: PGC-1 Blocking Peptide (301707)

PGC-1 Polyclonal Antibody (101707) PPARa Blocking Peptide (301710)

#### PPARδ LBD (human recombinant)

FAAR, NUC1, Nuclear Hormone Receptor 1, PPARB

**M.:** 54 kDa **Purity:** ≥95%

Supplied in: 50 mM sodium phosphate, pH 7.2, containing 20% glycerol, 150 mM sodium chloride, and 1 mN DTT

Source: Recombinant protein isolated from Baculovirus overexpression system in Sf21 cells

10 µg 25 µg

\*All 5-cis 2-series PGs (those containing a 5,6-double bond) will contain a small amount of the 5-trans isomer. This isomer is generally undetectable using normal phase silica columns and plates, but may be resolved using RP-HPLC. The purity for all such 2-series PGs excludes the 1-3% trans isomer which will generally be present.

PPAR Ligands					
Item No.	Product Name	Target	Mode of Action	Effective Concentration	Sizes
13452	AM3102	PPARα	Agonist	EC <sub>50</sub> = 100 nM	5 mg • 10 mg • 25 mg • 50 mg
60924	Azelaoyl PAF	PPARγ	Agonist	~ Equal to Rosiglitazone	500 μg • 1 mg • 5 mg • 10 mg
10009145	Bezafibrate	pan PPAR	Agonist	EC <sub>50</sub> = 20-60 μM	500 mg • 1 g • 5 g • 10 g
10009017	CAY10514	PPARα PPARγ	Dual Agonist	$EC_{50} = 0.173 \mu M (PPARα)$ $EC_{50} = 0.642 \mu M (PPARγ)$	1 mg • 5 mg • 10 mg • 50 mg
10008846	CAY10573	pan-PPAR	Agonist	$IC_{50} = 50-225 \mu\text{M}$	500 μg • 1 mg • 5 mg • 10 mg
10012536	CAY10592	PPARS	Agonist	EC <sub>50</sub> = 53 nM (rat)	1 mg • 5 mg • 10 mg • 50 mg
13282	CAY10599	PPARS	Agonist	EC <sub>50</sub> = 50 nM	1 mg • 5 mg • 10 mg • 25 mg
71730	Ciglitazone	PPARγ	Agonist	EC <sub>50</sub> = 3 μM	1 mg • 5 mg • 10 mg • 50 mg
10005745	Clofibrate	PPARα	Agonist	EC <sub>50</sub> = 55 μM (human)	500 μl • 1 ml • 5 ml • 10 ml
10005368	Fenofibrate	PPARα	Agonist	EC <sub>50</sub> = 30 μM (human)	1 g • 5 g • 10 g • 50 g
10006798	GW 0742	PPARS	Agonist	EC <sub>50</sub> = 1.1 nM	5 mg • 10 mg • 25 mg • 50 mg
10008613	GW 7647	PPARα	Agonist	EC <sub>50</sub> = 6 nM (human)	1 mg • 5 mg • 10 mg • 25 mg
10011211	GW 9578	PPARα	Agonist	EC <sub>50</sub> = 50 nM (human)	500 μg • 1 mg • 5 mg • 10 mg
70785	GW 9662	PPARy	Antagonist	Blocks differentiation of monocytes to osteoclasts by >90% at a dose of 0.1 $\mu M$	1 mg • 5 mg • 10 mg • 50 mg
10009880	GW 590735	PPARα	Agonist	$EC_{50} = 4 \text{ nM}$	1 mg • 5 mg • 10 mg • 25 mg
10009661	N-Octadecyl-N'-propyl-sulfamide	PPARα	Agonist	EC <sub>50</sub> = 100 nM	5 mg • 10 mg • 25 mg • 50 mg
90265	Oleoyl Ethanolamide	PPARα	Agonist	EC <sub>50</sub> = 120 nM	5 mg • 10 mg • 50 mg • 100 mg
18570	15-deoxy- $\Delta^{12,14}$ -Prostaglandin J $_2$	PPARS	Agonist	$EC_{50} = 7 \mu M$	100 μg • 500 μg • 1 mg • 5 mg
71740	Rosiglitazone	PPARγ	Agonist	$K_d = 43 \text{ nM}$	5 mg • 10 mg • 50 mg • 100 mg
10026	T0070907	PPARγ	Antagonist	$IC_{50} = 1$ nM for inhibition of Rosiglitazone binding	1 mg • 5 mg • 10 mg • 50 mg
71750	Troglitazone	PPARγ	Agonist	$EC_{50} = 0.55 \mu\text{M} \text{ (human)}$	5 mg • 10 mg • 50 mg • 100 mg
70730	Wy 14643	PPARα PPARδ	Agonist	Species dependent (PPARα: 0.1 - 10 μM)	5 mg • 10 mg • 50 mg • 250 mg

For a full product listing, please visit www.caymanchem.com

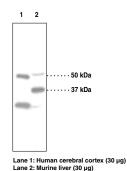
#### PPARδ Polyclonal Antibody

FAAR, NUC1, Nuclear Hormone Receptor 1, PPARβ

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

**Summary:** Antigen: human PPARδ amino acids 39-54 • Host: rabbit • Cross Reactivity: (+) human, murine, rat, ovine, and porcine PPARδ • Application(s): ICC, IHC, and WB • PPARδ is ubiquitously expressed but is particularly abundant in tissues such as liver, intestine, kidney, abdominal adipose, and skeletal muscle, all of which are involved in lipid metabolism. PPARδ is a mediator of diverse physiological functions including lipid and cholesterol homeostasis, embryo implantation, cancer development, and obesity.

500 µl



• Also Available: PGC-1 Blocking Peptide (301707)
PGC-1 Polyclonal Antibody (101707)
PPARδ Blocking Peptide (10006247)

PPARô Western Ready Control (10009568)

101720 PPARy FL (human recombinant from E. coli)

PPARy Full Length

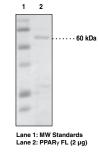
M<sub>r</sub>: ~60 kDa **Purity:** ≥90% by SDS-PAGE

Supplied in: 20 mM Tris HCl, pH 8.0, containing 250 mM potassium,

20% glycerol, 5 mM DTT, and 0.5 mM EDTA

Source: Human recombinant N-terminal His-tagged protein expressed in E. coli





61700

## PPAR<sub>Y</sub> FL (human recombinant from Sf21 cells)

10009987

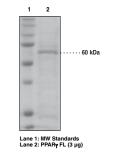
PPARy Full Length

M<sub>r</sub>: -60 kDa **Purity:** ≥80% by SDS-PAGE

**Supplied in:** 50 mM sodium phosphate, pH 7.2, containing 100 mM sodium chloride, 1 mM DTT, and 20% glycerol

Source: Human recombinant N-terminal His-tagged protein expressed in Sf21 insect cells





#### PPAR<sub>γ</sub> LBD (human recombinant)

10007941

PPARγ Ligand Binding Domain

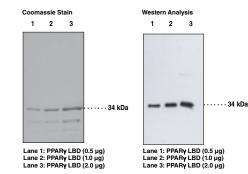
M<sub>r</sub>: ~34 kDa **Purity:** ≥90%

Supplied in:~50~mM~sodium phosphate, pH~7.2, containing~20%~glycerol, 150~mM

sodium chloride, and 1 mM DTT

 $\textbf{Source:} \ \text{Recombinant N-terminal His-tagged protein expressed in } \textit{E. coli}$ 





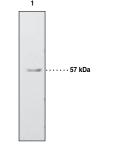
#### PPARy Polyclonal Antibody

101700

Peptide affinity-purified IgG **Stability:**  $\ge 1$  year at  $-20^{\circ}$ C

**Summary:** Antigen: human PPAR $\gamma_1$  amino acids 82-101 (amino acids 110-129 of PPAR $\gamma_2$ ) • Host: rabbit • Cross Reactivity: (+) human and murine PPAR $\gamma_1$  and PPAR $\gamma_2$  • Application(s): WB • PPAR $\gamma$  is a ligand-activated transcription factor involved in the regulation of lipid homeostasis and may function as a master regulator of adipogenesis.

500 μl



Lane 1: Rat adipose homogenate (30 µg)

\*Also Available: PGC-1 Blocking Peptide (301707)
PGC-1 Polyclonal Antibody (101707)
PPARγ Blocking Peptide (301700)

#### PPAR Transcription Factor Assay Kits

PPARs are ligand-activated transcription factors belonging to the large superfamily of nuclear receptors. PPAR $\alpha$  primarily activates genes encoding proteins involved in fatty acid oxidation, while PPAR $\gamma$  primarily activates genes directly involved in lipogenic pathway and insulin signaling. Members of the PPAR family are important direct targets of many antidiabetic and hypolipidemic drugs. Cayman's PPAR Transcription Factor Assays are a non-radioactive, sensitive method for detecting specific transcription factor DNA binding activity in nuclear extracts and whole cell lysates. A 96-well ELISA replaces the cumbersome radioactive electrophoretic mobility shift assay (EMSA). A specific dsDNA sequence containing the PPAR response element is immobilized onto the bottom of wells of a 96-well plate. PPARs contained in a nuclear extract, bind specifically to the PPAR response element. PPAR $\alpha$ ,  $\delta$ , or  $\gamma$  are detected by addition of specific primary antibodies directed against the individual PPARs. A secondary antibody conjugated to HRP is added to provide a sensitive colorimetric readout at 450 nm.

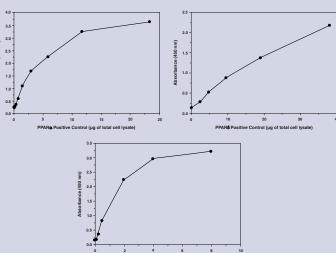
# PPAR $\alpha$ , $\delta$ , $\gamma$ Complete Transcription Factor Assay Kit

10008878

**Stability:** ≥1 year at -20°C

**Summary:**This kit contains individual primary antibodies for PPAR $\alpha$ ,  $\delta$ , or  $\gamma$  to follow detection of each receptor in separate wells of the plate.

96 wells

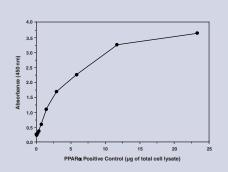


#### PPARα Transcription Factor Assay Kit

10006915

**Stability:** ≥6 months at -20°C

96 wells



• Also Available: PPARδ Transcription Factor Assay Kit (10006914)

PPARγ FP-Based Ligand Screening Assay Kit - Green (10007685)

PPARy Transcription Factor Assay Kit (10006855)

**36** Cayman Chemical caymanchem.com Cayman Chemical

# eNOS, Vascular Endothelial Health, and CVD



The vascular endothelium is an active organ. Its integrity supports the beneficial effects of antioxidant, anti-inflammatory, anticoagulant, and pro-fibrinolytic actions. Under normal circumstances, it releases nitric oxide (NO), which inhibits the adhesion and migration of leukocytes, the migration and proliferation of smooth muscle cells, the aggregation and adhesion of platelets, and is essential for maintaining vascular tone. Reduced NO activity leads to endothelial dysfunction, a systemic event that precedes the development of atherosclerosis and other cardiovascular complications (hypertension, congestive heart failure, stroke, etc.). Numerous studies have shown that functional endothelial NO synthase (eNOS), which synthesizes NO in the endothelium, is protective against these pathological vascular complications.<sup>1</sup> The synthesis of NO is a remarkably dynamic process of redox reactions and allosteric modifications, whereby a dysfunctional endothelium can be restored with appropriate therapeutics. Thus, understanding the signaling pathways that confer vascular protection is of central importance to the refinement of therapeutic approaches to prevent or ameliorate cardiovascular diseases. What follows is a brief overview of eNOS signaling and protein modifications critical to proper endothelial function.

#### **eNOS Activation and NO Signaling**

In endothelial cells, functional eNOS transfers electrons from NADPH, via the flavins FAD and FMN in the carboxy-terminal reductase domain, to the heme in the amino-terminal oxygenase domain. (Figure 1) Here the substrate L-arginine is oxidized to L-citrulline and NO. eNOS is indirectly activated by physiological and metabolic stimuli, laminar shear stress, and receptor-dependent agonists (i.e., VEGF, PAF, bradykinin). eNOS is bound to caveolin-1 through a consensus site in caveolae, which regulates the basal state of NO production. Its activation involves dissociation from caveolin-1, association with heat shock protein (Hsp)-90, Ca<sup>2+</sup>- calmodulin binding, (calmodulin-dependent kinase II, Akt, PKA, or AMPK) phosphorylation of serine-1177, and dephosphorylation of threonine-495. At least six phosphorylation sites have been indentified for roles in eNOS activity. These include serine-116, -617, -635, and -1179 as well as threonine-497 and tyrosine-83. Calcium-calmodulin activation and phosphorylation increase NO production.

In endothelial cells eNOS is found mainly in the plasma membrane and the Golgi complex, but it also distributes in the cytosol. eNOS bound to the plasma membrane releases greater amounts of basal NO compared to that with intracellular localization, resulting in a greater ability to elicit cGMP signaling.2 However, translocation of eNOS to subcellular compartments may be useful for precisely directing an efficacious signal to its intended target before encountering scavengers of reactive species. Movement from plasma membrane to subcellular locations (including the Golgi) is associated with de-myristoylation and de-palmitoylation, which encourage dissociation of eNOS with caveolin-1. Caveolae either pinch off (endocytosis) the plasma membrane, internalizing eNOS into the cell, or dissociate in some other manner.

Once NO is produced by the endothelium, it can regulate various aspects of the vascular system by targeting soluble guanylyl cyclase (sGC), which mediates cGMP-PKG signaling, or it can initate nitrosation reactions with iron-sulphur-centered proteins or proteins with reactive thiols (S-nitrosylation).(Figure 2) The PKG-dependent signaling pathway induces relaxation of vascular smooth muscle and suppression of platelet aggregation. Protein S-nitrosylation, a post-translational modification that has been compared to phosphorylation where NO modifies cysteine residues of target proteins, induces allosteric changes in protein function. It participates in the regulation of apoptosis, proliferation, protein secretion, ion channel activity, vesicular trafficking of local proteins, and NOS activity. Location of eNOS can influence these signaling events. eNOS activity in the Golgi results in the compartmentalization of NO and S-nitrosylation of select local proteins. However, Golgi-restricted eNOS has also been demonstrated to supply NO to adjacent smooth muscle cells, contributing to endothelium-dependent relaxation.<sup>2</sup> The factors are not yet clear as to what distinguishes signaling through cGMPdependent vs. S-nitrosylation modifications.

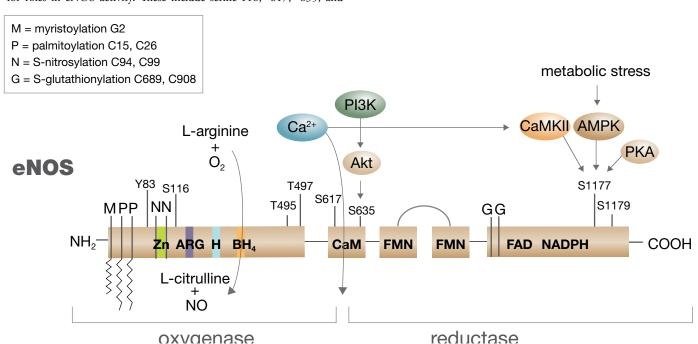


Figure 1. eNOS Protein Structure. Key phosphorylation sites are numbered and additional modifications are indicated by an M (myristoylation), P (palmitoylation), N (S-nitrosylation), or G (S-glutathionylation).

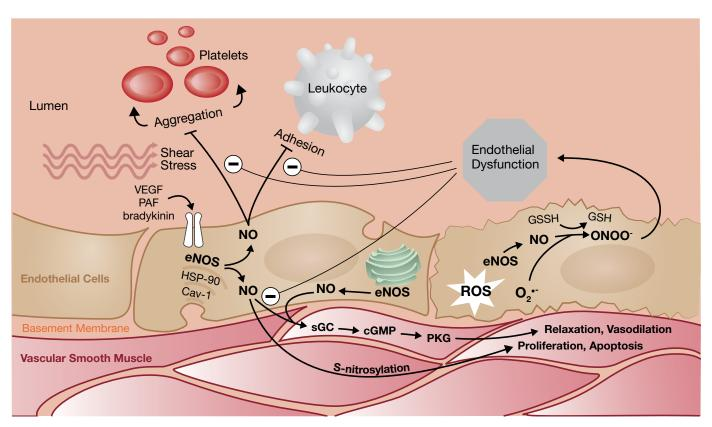


Figure 2. Schematic Representation of Endothelial Nitric Oxide Signaling. eNOS is activated by physiological and metabolic stimuli, shear stress, and receptor-dependent agonists to functionally inhibit platelet aggregation, leukocyte adhesion, and smooth muscle cell proliferation, as well as to maintain vascular tone. Oxidative stress uncouples eNOS-derived

#### **Uncoupling of eNOS Function**

Cardiovascular risk factors induce oxidative stress, which alters various functions of the endothelium, including the reduced bioavailability of NO and the increased production of superoxide (O<sub>2</sub>·-) and reactive oxygen species (ROS) in the vessel wall. These risk factors lead to an enhanced production of NADPH oxidases. NADPH-oxidase-derived superoxide readily reacts with eNOS-derived NO to form peroxynitrite (ONOO<sup>-</sup>). The essential eNOS cofactor (6R-)5,6,7,8-tetrahydrobiopterin (BH<sub>4</sub>) is highly sensitive to oxidation by ONOO-. With BH<sub>4</sub> deficiency, reduction of O<sub>2</sub> becomes uncoupled from NO synthesis, thereby converting eNOS to a superoxide-producing enzyme.

Recently, the uncoupling of oxygen reduction by eNOS (i.e., the switch from NO synthesis to superoxide production) has been proposed to result from S-glutathionylation of eNOS. This is a post-translational modification of key cysteine residues where glutathione is added through reversible thiol-disulfide exchange with oxidized glutathione (GSSH) or reaction of oxidant-induced protein thiyl radicals with reduced glutathione (GSH). S-glutathionylation of eNOS at cysteine-689 and cysteine-908 reversibly decreases NOS activity while at the same time increases superoxide generation.<sup>3,4</sup> S-glutathionylation of eNOS is increased with impaired endothelium-dependent vasodilation in hypertensive vessels.<sup>3,4</sup> Restoration of this function has been achieved in an in vitro model with thio-specific reducing agents that reverse S-glutathionylation of eNOS.<sup>3,4</sup>

#### Therapeutic Potential

Compounds interfering with the renin-angiotensin-aldosterone system as well as statins can reduce vascular oxidative stress and increase bioactive NO. Additionally, new small molecules are being identified that have the potential to prevent eNOS uncoupling and, at the same time, enhance eNOS expression. Compounds that increase eNOS protein levels are only advantageous if they support eNOS functionality otherwise superoxide production would be reinforced. The protein kinase C inhibitor PKC 412 (also known as midostaurin; Item No. 10459) maintains eNOS functionality in disease while at the same time upregulating the expression of the enzyme. It does so though a PKC-independent mechanism and reduces NADPH oxidase expression via PKC inhibition. By reducing NADPH oxidase-mediated oxidative stress, PKC 412 reverses eNOS uncoupling in spontaneously hypertensive rats and in atherosclerosisprone apolipoprotein E knockout mice.<sup>5,6</sup> The polyphenolic phytoalexin trans-resveratrol (Item No. 70675) also reverses eNOS uncoupling and enhances eNOS expression. It has been shown to upregulate superoxide dismutases, glutathione peroxidase 1, and catalase in hypercholesterolemic apolipoprotein E knockout mice. Additionally it downregulates NADPH oxidases resulting in an overall reduction of peroxynitrite levels.<sup>5,7,8</sup> The continued development of compounds that combine eNOS upregulation with eNOS recoupling may thus have therapeutic potential for cardiovascular diseases and the restoration of proper endothelial function.

- Atochin, D.N. and Huang, P.L. Eur. J. Physiol. 460, 965-974 (2010).
- Qian, J., Zhang, Q., Church, J.E., et al. Am. J. Physiol. Heart Circ. Physiol. 298, H112-H118 (2010).
- Duan, D.D. and Kwan, C-Y. Acta Pharmacol. Sin. [Epub ahead of print] (2011) Chen, C-A., Wang, T-Y., Varadharaj, S., et al. Nature 468, 1115-1120 (2010)
- Li, H., Hergert, S.M., Schäfer, S.C., et al. Nitric Oxide 12, 231-236 (2005).
- Li. H. and Förstermann. U. Meeting contribution at Deutsche Gesellschaft für Arterioskleroseforschung. Blaubeuren
- Xia, H., Daiber, A., Habermeier, A., et al. J. Pharmacol. Exp. Ther. 335(1), 149-154 (2010).
- Spanier, G., Xu, H., Xia, N. et al. J. Physiol. Pharmacol. 60(Suppl 4), 111-116 (2009)

#### Pravastatin (sodium salt)

10010343

10007599

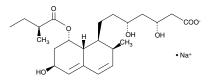
515121

[81131-70-6]

MF: C<sub>23</sub>H<sub>35</sub>O<sub>7</sub> • Na FW: 446.5 Purity: ≥98% A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent, competitive inhibitor of HMG-CoA reductase (K<sub>i</sub> = 2.3 nM)

10 mg 25 mg 50 mg



#### Prorenin (human recombinant)

M.: 45 kDa Purity: ≥99% by SDS-PAGE **Supplied in:** 50 mM mM Tris, pH 8.0

Source: Recombinant enzyme expressed in HEK cells

50 µg 100 µg 500 µa

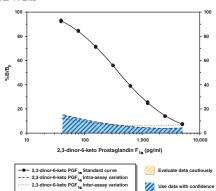
#### 2,3-dinor-6-keto Prostaglandin F<sub>1a</sub> EIA Kit

**Stability:** ≥1 year at -20°C

Sensitivity: 50% B/B<sub>0</sub>: 400 pg/ml • 80% B/B<sub>0</sub>: 100 pg/ml

Summary: PGI, is non-enzymatically hydrated to 6-keto  $PGF_{1\alpha}$ , and then quickly converted to the major urinary metabolite, 2,3-dinor-6-keto PGF<sub>10</sub>. The majority of 6-keto PGF<sub>10</sub> in urine is of renal origin with only 14% originating from plasma. Cayman's 2,3-dinor-6-keto PGF<sub>10</sub> EIA utilizes a highly selective monoclonal antibody that exhibits no cross reactivity with 6-keto PGF<sub>10</sub>, thus providing a method for accurate measurement of systemic PGI<sub>2</sub> production.

96 strip/solid wells 480 strip/solid wells



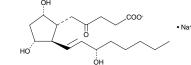
#### 2,3-dinor-6-keto Prostaglandin $F_{1\alpha}$ (sodium salt) 15120

MF: C<sub>10</sub>H<sub>20</sub>O<sub>6</sub> • Na FW: 364.4 Purity: ≥98%

A solution in methanol **Stability:** ≥2 years at -20°C

**Summary:** β-oxidation product of 6-keto PGF<sub>1α</sub> and the major urinary metabolite of PGI<sub>2</sub> in humans





• Also Available: 2,3-dinor-6-keto Prostaglandin F<sub>10</sub>-d<sub>9</sub> (sodium salt) (9000462)

6-keto Prostaglandin F<sub>1</sub>

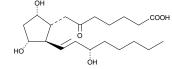
[58962-34-8]

**MF:**  $C_{20}H_{34}O_6$  **FW:** 370.5 **Purity:**  $\geq$ 99%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: Inactive, non-enzymatic hydrolysis product of PGI2; serves as a useful marker of PGI2 biosynthesis in vivo

5 mg 10 mg 50 mg



• Also Available: 6-keto Prostaglandin F<sub>10</sub>-d<sub>4</sub> (315210)

6-keto Prostaglandin F<sub>10</sub> Lipid Maps MS Standard (10007219)

6-keto Prostaglandin  $F_{1\alpha}$ -d<sub>4</sub> Lipid Maps MS Standard (10007274)

6-keto Prostaglandin F<sub>10</sub> Quant-PAK (10006830)

#### 6-keto Prostaglandin F<sub>1</sub>, EIA Kit

515211

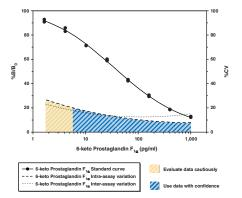
15210

**Stability:** ≥1 year at -20°C

**Sensitivity:** 50% B/B<sub>0</sub>: 40 pg/ml • 80% B/B<sub>0</sub>: 6 pg/ml

Summary: Prostacyclin is formed from arachidonic acid primarily by the vascular endothelium and renal cortex. It is a potent vasodilator and inhibitor of platelet aggregation. PGI<sub>2</sub> is non-enzymatically hydrated to 6-keto PGF<sub>10</sub> ( $t_{1/2} = 2-3$ minutes), and then quickly converted to the major metabolite, 2,3-dinor-6-keto  $PGF_{1\alpha}$  ( $t_{1/2} = 30$  minutes).

96 strip/solid wells 480 strip/solid wells



#### 9,11-methane-epoxy Prostaglandin F<sub>1</sub>,

10007850

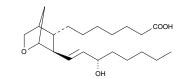
[72517-81-8] 9,11-epoxymethano PGH<sub>1</sub>

MF:  $C_{21}H_{36}O_4$  FW: 352.5 Purity:  $\geq 96\%$ 

A solution in methyl acetate **Stability:** ≥1 year at -20°C

Summary: A stable epoxymethano analog of PGH<sub>1</sub>; produces a strong and doserelated aggregation of washed rabbit platelets (EC<sub>50</sub> =  $0.88 \mu M$ )

100 µg 500 µg 1 mg 5 mg



#### Prostaglandin I Synthase Monoclonal Antibody (Clone isn-1)

160630

PGIS, PGI Synthase, Prostacyclin Synthase

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

**Summary:** Antigen: bovine lung PGIS • Host: mouse, clone isn-1 • Cross Reactivity: (+) bovine, murine, rat, ovine, guinea pig, and rabbit PGIS • Isotype: IgG<sub>1</sub> • Application(s): IHC and IP; WB not recommended • PGIS catalyzes the conversion of PGH2 to PGI2 (prostacyclin).

1 ea

#### Prostaglandin I Synthase Polyclonal Antibody 160640

PGIS, PGI Synthase, Prostacyclin Synthase

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: bovine PGIS amino acids 299-329 • Host: rabbit • Cross Reactivity: (+) bovine, ovine, and human PGIS; (-) rat PGIS • Application(s): IP and WB • PGIS catalyzes the isomerization of PGH2 to PGI2, a potent vasodilator and inhibitor platelet aggregation.

500 µl

• Also Available: Prostaglandin I Synthase Blocking Peptide (360640)

#### Prostaglandin I Synthase (murine) Polyclonal Antibody

100023

18220

PGIS, PGI Synthase, Prostacyclin Synthase

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: murine PGIS amino acids 475-490 conjugated to KLH • Host: rabbit • Cross Reactivity: (+) human, bovine, ovine, rat, and murine PGIS • Application(s): IHC and WB • PGIS catalyzes the isomerization of PGH2 to PGI2 (prostacyclin), a potent vasodilator and inhibitor of platelet aggregation.

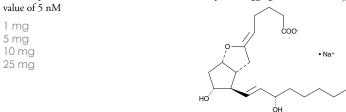
• Also Available: Prostaglandin I Synthase (murine) Blocking Peptide (10005257)

#### Prostaglandin I<sub>2</sub> (sodium salt)

[61849-14-7] Epoprostenol (sodium salt), Prostacyclin (sodium salt)

MF:  $C_{20}H_{31}O_5 \bullet \text{ Na FW: } 374.5 \text{ Purity: } \ge 99\%^*$ A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A primary metabolite of arachidonic acid formed by COX and PGIS; acts as a potent vasodilator and inhibitor of human platelet aggregation with an IC<sub>50</sub>



• Also Available: Prostaglandin I<sub>3</sub> (sodium salt) (18300)

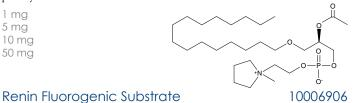
#### Pyrrolidino PAF C-16 60909

[91021-63-5]

MF: C<sub>28</sub>H<sub>56</sub>NO<sub>7</sub>P FW: 549.7 Purity: ≥98%

A lyophilized powder **Stability:** ≥1 year at -20°C

Summary: A synthetic analog of PAF C-16 with approximately 3-10 times increased potency



#### **Purity:** ≥95%

A red/brown crystalline solid **Stability:** ≥1 year at -20°C

**Summary:** The renin fluorogenic substrate consists of the normal peptide substrate for renin which has been linked to the fluorophore EDANS at one end and to a nonfluorescent quenching molecule (Dabcyl) at the other. After cleavage by renin, the product (peptide-EDANS) is brightly fluorescent and can be easily analyzed using an excitation wavelength of 340 nm and emission wavelengths of 485-510 nm.

1 mg

#### 10006217 Renin (human recombinant)

M.: -40 kDa Purity: ≥99% by SDS-PAGE

**Supplied in:** Sodium acetate buffer **Stability:** ≥6 months at -80°C

**Source:** Recombinant enzyme expressed in HEK cells as prorenin. The prorenin was activated using trypsin and purified using peptide affinity chromatography.

10 µg

25 µg

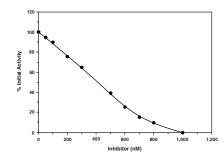
50 µg

Renin Inhibitor Screening Assay Kit

10006270

**Stability:** ≥6 months at -80°C

Summary: Cayman's Renin Inhibitor Screening Assay provides a convenient 96-well format for evaluating human renin inhibitors. The assay utilizes a renin-based synthetic peptide substrate which incorporates the fluorophore EDANS at one end and an EDAN-quenching molecule (Dabcyl) at the other end. After cleavage by renin, the peptide-EDANS product is released yielding bright fluorescence.



#### Resorcinonaphthalein

13082

[41307-63-5] NSC 354317

**MF:**  $C_{24}H_{14}O_5$  **FW:** 382.4 **Purity:**  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: Increases the activity of angiotensin-converting enzyme 2 (ACE2) in vitro  $(EC_{50} = 19.5 \mu M)$ 

5 mg 10 mg

50 mg 100 mg

#### Ro 48-8071 10006415

[161582-11-2]

**MF:** C<sub>23</sub>H<sub>27</sub>BrFNO<sub>2</sub> **FW:** 448.4 **Purity:** ≥98%

A solution in methyl acetate **Stability:** ≥1 year at -20°C

Summary: An inhibitor of oxidosqualene cyclase (OSC) that has LDL cholesterol lowering activity; inhibits OSC from human liver microsomes and HepG2 cells with IC<sub>50</sub> values of approximately 6.5 and 1.5 nM, respectively

5 mg 10 mg 50 mg

<sup>\*</sup>All 5-cis 2-series PGs (those containing a 5,6-double bond) will contain a small amount of the 5-trans isomer. This isomer is generally undetectable using normal phase silica columns and plates, but may be resolved using RP-HPLC. The purity for all such 2-series PGs excludes the 1-3% trans isomer which will generally be present.

#### 10010344 Simvastatin

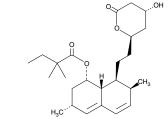
[79902-63-9] MK 733

**MF:** C<sub>25</sub>H<sub>38</sub>O<sub>5</sub> **FW:** 418.6 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** A potent, competitive inhibitor of HMG-CoA reductase ( $K_i = 0.12 \text{ nM}$ )

10 mg 25 mg 50 mg



	Epoxy Fatty Acids				
Item No.	Product Name	Sizes			
Arachido	nic Acid Derivatives				
50211	(±)5(6)-EET	25 μg • 50 μg • 100 μg • 500 μg			
10009984	(±)5(6)-EET-d <sub>11</sub>	10 µg • 25 µg • 50 µg • 100 µg			
10008596	(±)5(6)-EET Ethanolamide	25 μg • 50 μg • 100 μg • 250 μg			
50210	(±)5(6)-EET methyl ester	25 µg • 50 µg • 100 µg • 250 µg			
50351	(±)8(9)-EET	25 µg • 50 µg • 100 µg • 500 µg			
10009532	(±)8(9)-EET-d <sub>11</sub>	10 µg • 25 µg • 50 µg • 100 µg			
10008597	(±)8(9)-EET Ethanolamide	25 µg • 50 µg • 100 µg • 250 µg			
50350	(±)8(9)-EET methyl ester	25 µg • 50 µg • 100 µg • 250 µg			
50511	(±)11(12)-EET	25 μg • 50 μg • 100 μg • 500 μg			
10006413	(±)11(12)-EET-d <sub>11</sub>	10 µg • 25 µg • 50 µg • 100 µg			
10008598	(±)11(12)-EET Ethanolamide	25 μg • 50 μg • 100 μg • 500 μg			
50510	(±)11(12)-EET methyl ester	25 μg • 50 μg • 100 μg • 500 μg			
50651	(±)14(15)-EET	25 μg • 50 μg • 100 μg • 500 μg			
10006410	(±)14(15)-EET-d <sub>11</sub>	25 µg • 50 µg • 100 µg • 250 µg			
10008599	(±)14(15)-EET Ethanolamide	25 µg • 50 µg • 100 µg • 250 µg			
50650	(±)14(15)-EET methyl ester	25 μg • 50 μg • 100 μg • 250 μg			
10009286	(±)14(15)-EET-SI	25 μg • 50 μg • 100 μg • 250 μg			
Docosahe	exanenoic Acid Derivatives				
90314	4(5)-EpDPE methyl ester	25 µg • 50 µg • 100 µg • 500 µg			
10174	16(17)-EpDPE	25 μg • 50 μg • 100 μg • 500 μg			
10175	19(20)-EpDPE	25 µg • 50 µg • 100 µg • 500 µg			
Eicosape	ntaenoic Acid Derivatives				
90114	5(6)-EpETE methyl ester	25 µg • 50 µg • 100 µg • 500 µg			
10173	14(15)-EpETE	25 μg • 50 μg • 100 μg • 500 μg			
50861	17(18)-EpETE	25 µg • 50 µg • 100 µg • 500 µg			
Linoleic A	Acid Derivatives				
52400	(±)9(10)-Ep0ME	25 μg • 50 μg • 100 μg • 250 μg			
10009995	(±)9(10)-Ep0ME-d <sub>4</sub>	25 µg • 50 µg • 100 µg • 250 µg			
52450	(±)12(13)-Ep0ME	25 µg • 50 µg • 100 µg • 250 µg			
10009996	(±)12(13)-Ep0ME-d <sub>4</sub>	25 µg • 50 µg • 100 µg • 250 µg			

Also Available:  $(\pm)$ 5(6)-EET Lipid Maps MS Standard (10007260) •  $(\pm)$ 8(9)-EET Lipid Maps MS Standard (10007261) • (±)11(12)-EET Lipid Maps MS Standard (10007262) •

(±)14(15)-EET Lipid Maps MS Standard (10007263)

#### Soluble Epoxide Hydrolase Cell-Based Assay Kit

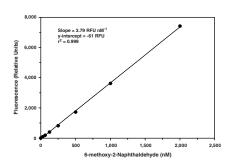
600090

CEH, Cytosolic Epoxide Hydrolase, EPHX2, Epoxide Hydrolase 2

**Stability:** ≥1 year at -80°C

**Summary:** Mammalian sEH is a member of the  $\alpha/\beta$ -hydrolase fold family of enzymes that catalyze the hydrolysis of exogenous and endogenous epoxides to vicinal diols. Endogenous substrates for sEH include EETs which exhibit vasodilatory and antiinflammatory activity. Cayman's sEH Cell-Based Assay provides a fluorescence-based method for detecting epoxide hydrolase activity in whole cells. The assay utilizes Epoxy Fluor 7, a sensitive fluorescent substrate for sEH that can be used to monitor the activity of both human and murine enzymes.

480 wells



#### Soluble Epoxide Hydrolase (human recombinant)

10011669

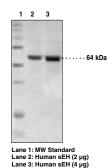
CEH, Cytosolic Epoxide Hydrolase, EPHX2, Epoxide Hydrolase 2

**M**<sub>r</sub>: 67 kDa **Purity**: ≥95%

Supplied in: TBS buffer at pH 7.4 containing 20% glycerol

**Source:** Recombinant N-terminal His-tagged protein purified from Sf21 cells

50 µg 100 µg



#### Soluble Epoxide Hydrolase Inhibitor Screening Assay Kit

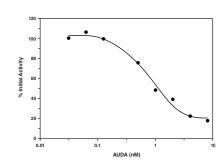
10011671

CEH, Cytosolic Epoxide Hydrolase, EPHX2, Epoxide Hydrolase 2

**Stability:** ≥6 months at -80°C

Summary: sEH catalyzes the hydrolysis of exogenous and endogenous epoxides to vicinal diols. Endogenous substrates for sEH include epoxyeicosatrienoic acids (EETs) which exhibit vasodialatory and anti-inflammatory activity. Small molecule inhibitors of sEH may therefore hold promise as therapeutics for the treatment of hypertension and vascular inflammation. Cayman's fluorescence-based sEH Inhibitor Screening Assay provides a convenient method for screening epoxide hydrolase inhibitors. The assay utilizes (3-phenyl-oxiranyl)-acetic acid cyano-(6-methoxy-naphthalen-2-yl)methyl ester (PHOME) as a substrate. Hydrolysis of PHOME by epoxide hydrolase produces the highly fluorescent 6-methoxy-2-naphthaldehyde which can be analyzed using an excitation wavelength of 330 nm and emission wavelength of 465 nm. Each kit contains buffer, substrate, and sufficient human recombinant sEH for 100 tests.

96 wells



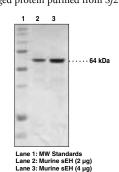
#### Soluble Epoxide Hydrolase (murine recombinant from Sf21 cells)

10011670

CEH, Cytosolic Epoxide Hydrolase, EPHX2, Epoxide Hydrolase 2 M<sub>r</sub>: 64 kDa Purity: ≥95%

Supplied in: TBS buffer at pH 7.4 containing 20% glycerol **Source:** Recombinant N-terminal His-tagged protein purified from Sf21 cells

50 µg 100 µg

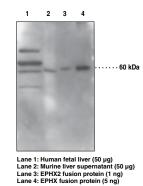


#### Soluble Epoxide Hydrolase Polyclonal Antibody

10010146

CEH, Cytosolic Epoxide Hydrolase, EPHX2, Epoxide Hydrolase 2 Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

**Summary:** Antigen: rat sEH amino acids 292-306 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat sEH • Application: WB • sEH converts epoxyeicosatrienoic acids (EpETrEs, EETs) to the corresponding dihydroxy eicosatrienoic acids (DiHETrEs, DHETs), which diminishes their vasodilator activity.



• Also Available: Soluble Epoxide Hydrolase Blocking Peptide (10010147)

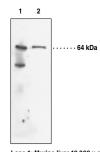
#### Soluble Epoxide Hydrolase (FL) Polyclonal Antibody

13560

CEH, Cytosolic Epoxide Hydrolase, EPHX2, Epoxide Hydrolase 2 Protein A-purified IgG **Stability:** ≥1 year at -20°C

**Summary:** Antigen: murine recombinant sEH • Host: rabbit • Cross Reactivity: (+) human and murine sEH • Application(s): WB • sEH converts epoxyeicosatrienoic acids (EpETrEs, EETs) to the corresponding dihydroxy eicosatrienoic acids (DiHETrEs, DHETs), which diminishes their vasodilator activity.

500 µl



supernatant Lane 2: Purified murine sEH

Dihydroxy Eicosatetraenoic Acids (DiHETEs)			
Item No.	Product Name	Sizes	
Arachidon	ic Acid Derivatives		
35200	5(S),6(R)-DiHETE	25 µg • 50 µg • 100 µg • 250 µg	
10009628	5(S),6(R)-11- <i>trans</i> -DiHETE	25 µg • 50 µg • 100 µg • 250 µg	
35210	5(S),6(S)-DiHETE	25 µg • 50 µg • 100 µg • 250 µg	
35280	5(S),15(S)-DiHETE	25 µg • 50 µg • 100 µg • 250 µg	
35370	8(S),15(S)-DiHETE	25 µg • 50 µg • 100 µg • 250 µg	
Eicosapen	taenoic Acid Derivatives		
10006998	(±)14,15-DiHETE	25 µg • 50 µg • 100 µg • 250 µg	

10006999	(±)17,18-DiHETE	25 μg • 50 μg • 100 μg • 250 μg
Also Availa	able: 5(S),6(R)-DiHETE Lipid Maps M	S Standard (10007252) • 5(S),6(S)-DiHET
Lipid Ma	os MS Standard (10007253) • 5(S),1	5(S)-DiHETE Lipid Maps MS Standard

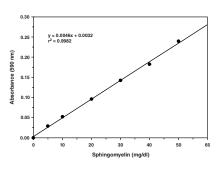
(10007255)

Sphingomyelin Assay Kit 10009928

**Stability:** ≥6 months at 4°C

Summary: Sphingomyelin is an important lipid component of cell membranes and lipoproteins. Cayman's SM Assay provides a specific, sensitive, and convenient method for quantifying SM in plasma or serum in a 96-well format with a colorimetric readout at 595 nm.

96 wells



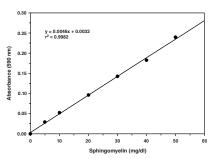
#### Sphingomyelinase Assay Kit

10006964

**Stability:** ≥6 months at -20°C

Summary: Cayman's Sphingomyelinase Assay provides a simple, reproducible, and sensitive tool for measuring neutral and acidic sphingomyelinase activity from tissue homogenates, cell lysates, serum, saliva, and urine. The assay employes a coupled enzymatic reaction system, resulting in the formation of highly fluorescent resorufin (excitation: 530-540 nm; emission 585-595 nm).

96 wells



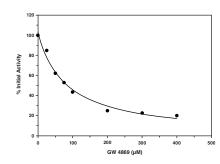
#### Sphingomyelinase Inhibitor Screening Assay Kit

700330

**Stability:** ≥6 months at -20°C

Summary: Cayman's Sphingomyelinase Inhibitor Screening Assay is a novel method for screening sphingomyelinase inhibitors. Cleavage of a unique sphingomyelin conjugate by SMase in the assay results in the release of a ceramide analog containing a free thiol which is detected by the SMase Detector to yield a highly fluorescent product.

96 wells



SQ 29,548 19025

[98672-91-4]

**MF:**  $C_{21}H_{29}N_3O_4$  **FW:** 387.5 **Purity:**  $\geq$ 99%\*

A crystalline solid **Stability:** ≥1 year at -20°C

**Summary:** A highly selective TP receptor antagonist  $(K_i = 4.1 \text{ nM})$ 

5 mg 10 mg 100 mg

#### SREBP-2 Cell-Based Translocation Assay Kit

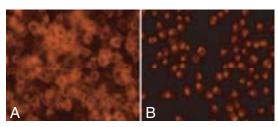
10009239

SREBF-2

**Stability:** ≥1 year at -20°C

Summary: Cayman's SREBP-2 Cell-Based Translocation Assay Kit provides the tools needed to study SREBP-2 movement within whole cells. The kit contains a highly specific SREBP-2 primary antibody together with a DyLight™ (trademarked by Pierce Biotechnology Inc.) conjugated secondary antibody in a ready to use format. Also included as a positive control is a cholesterol trafficking inhibitor, U18666A, which has been shown to activate SREBP-2 translocation into nuclei by scientists at Cayman Chemical Company.

96 wells



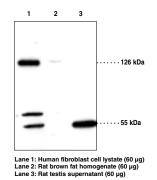
Translocation of SREBP-2 into nuclei by 24 µM U-18666A. Raw 264.7 cells were treated with DMSO (vehicle) or 24  $\mu$ M U-18666A for 72 hours. Panel A: Cells treated with DMSO alone demonstrate cytoplasmic localization of SREBP-2, indicating that most of cells have inactive protein. Panel B: U-18666A treatment for three days induced SREBP-2 translocation into the nuclei, indicating that blockage of cholesterol transport in these cells activates the protein.

#### SREBP-2 Polyclonal Antibody

10007663

Affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human SREBP-2 amino acids 455-469 • Host: rabbit • Cross Reactivity: (+) human, murine, and rat SREBP-2 • Application(s): ICC and WB • SREBP-2 is a transcription factor that plays a critical role in lipid homeostasis by regulating genes involved in cholesterol and fatty acid metabolism.



• Also Available: SREBP-2 Blocking Peptide (10009266) SREBP-2 Western Ready Control (10009749)

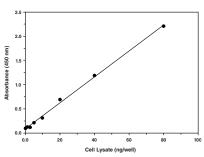
#### SREBP-2 Transcription Factor Assay Kit

**Stability:** ≥6 months at -20°C

Summary: SREBP-2 is a transcription factor that performs a critical role in the transcriptional regulation of genes involved in cholesterol synthesis and uptake including HMG-CoA synthase, HMG-CoA reductase, and the LDL receptor. Cayman's SREBP-2 Transcription Factor Assay is a non-radioactive, sensitive method for detecting SREBP-2 DNA binding activity in nuclear extracts and whole cell

96 wells

SREBF-2



T0901317 71810

[293754-55-9]

MF:  $C_{17}H_{13}F_0NO_2S$  FW: 481.3 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥1 year at -20°C

Summary: A potent and selective agonist for both LXRα and LXRβ (EC<sub>50</sub> =

5 mg 10 mg 50 mg 100 ma

10011348

#### Taprostene (sodium salt)

[87440-45-7] CG 4203, Rheocyclan

**MF:**  $C_{24}H_{29}O_5$  • Na **FW:** 420.5 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A stable prostacyclin analog and agonist of the IP receptor; does not activate EP4

500 µg 1 mg 5 mg 10 ma

# 90500

#### 3-Thiatetradecanoic Acid

[116296-31-2] 3-TDA

MF:  $C_{13}H_{26}O_2S$  FW: 246.4 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An analog of myristic acid wherein the C-3 carbon has been replaced by sulfur in a thioether linkage; acts as a PPAR ligand, increases fatty acid oxidation, and lowers plasma lipid levels

1 mg 5 mg 10 mg 50 ma

#### Thromboxane B<sub>2</sub>

19030

[54397-85-2]

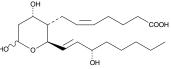
10007819

MF:  $C_{20}H_{34}O_6$  FW: 370.5 Purity:  $\geq 99\%$ \*

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A stable, biologically inert metabolite formed from the rapid nonenzymatic hydrolysis of TXA2

5 mg 10 mg 50 mg



• Also Available: Thromboxane B<sub>2</sub>-d<sub>4</sub> (319030)

Thromboxane B<sub>2</sub> Lipid Maps MS Standard (10007237) Thromboxane B<sub>2</sub> Quant-PAK (10006832)

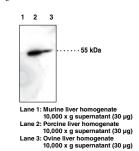
#### Thromboxane B<sub>2</sub> 11-dehydrogenase Polyclonal Antiserum

160720

Lyophilized antiserum **Stability:** ≥3 years at -20°C

Summary: Antigen: human erythrocyte TXB<sub>2</sub> 11-dehydrogenase • Host: rabbit • Cross Reactivity: (+) human erythrocyte, porcine liver, and ovine liver TXB2 11-dehydrogenase • Application(s): WB • TXB<sub>2</sub> 11-dehydrogenase catalyzes the conversion of TXB<sub>2</sub> to 11-dehydro TXB<sub>2</sub>

500 µl



#### Thromboxane B<sub>2</sub> EIA Kit

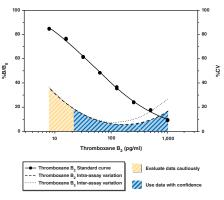
519031

**Stability:** ≥1 year at -20°C

**Sensitivity:** 50% B/B<sub>0</sub>: 57 pg/ml • 80% B/B<sub>0</sub>: 11 pg/ml

**Summary:** TXA<sub>2</sub> is produced from arachidonic acid by many cells and causes irreversible platelet aggregation and vascular and bronchial smooth muscle contraction. TXA2 is rapidly hydrolyzed non-enzymatically to form TXB2, which is then quickly metabolized ( $t_{1/2} = 5-7$  minutes) to urinary metabolites for clearance by the kidneys. Because of the transient nature of this compound it is difficult to accurately measure circulating levels in whole-animal experimental models. In fact, it has been shown that plasma and urine levels of TXB2 are primarily due to ex vivo platelet activation and intra-renal production, respectively. Therefore, measurement of TXB<sub>2</sub> metabolites such as 11-dehydro TXB<sub>2</sub> (Item No. 519501) and 2,3-dinor TXB2 (Item No. 519051) in urine and plasma may give better estimates of in vivo TXA<sub>2</sub> production. TXB<sub>2</sub> measurement is better suited towards samples that are not expected to undergo extensive metabolism such as perfusates, lavage samples, and tissue/cell culture medium or lysates.

96 strip/solid wells 480 strip/solid wells



• Also Available: Thromboxane B2 Express EIA Kit - Monoclonal (10004023)

<sup>\*</sup>All 5-cis 2-series PGs (those containing a 5,6-double bond) will contain a small amount of the 5-trans isomer This isomer is generally undetectable using normal phase silica columns and plates, but may be resolved using RP-HPLC. The purity for all such 2-series PGs excludes the 1-3% trans isomer which will generally be present.

#### 11-dehydro Thromboxane B<sub>2</sub>

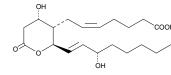
[67910-12-7] 11-keto TXB<sub>2</sub>

MF:  $C_{20}H_{32}O_6$  FW: 368.5 Purity:  $\geq 98\%^*$ 

A solution in methyl acetate **Stability:** ≥1 year at -20°C

Summary: One of the main plasma and urinary metabolites of TXB2 that is used as a marker of TXA2 synthesis in vivo

100 µg 250 µg 500 µg 1 mg



• Also Available: 11-dehydro Thromboxane B<sub>2</sub>-d<sub>4</sub> (319500)

11-dehydro Thromboxane B, Lipid Maps MS Standard (10007239)

11-dehydro Thromboxane B, Quant-PAK (10006832)

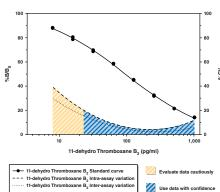
#### 11-dehydro Thromboxane B2 EIA Kit

**Stability:** ≥1 year at -20°C

Sensitivity: 50% B/B<sub>0</sub>: 93 pg/ml • 80% B/B<sub>0</sub>: 16 pg/ml

Summary: TXA2 is produced from arachidonic acid by many cells and causes irreversible platelet aggregation and vascular and bronchial smooth muscle contraction. TXA2 is rapidly hydrolyzed non-enzymatically to form TXB2. Although it is common to estimate TXA2 levels by measuring TXB2, most of the TXB<sub>2</sub> measured in plasma or urine is due to ex vivo platelet activation or intrarenal production, respectively. Measurement errors are compounded by the fact that normal concentrations of circulating TXB, are extremely low (1-2 pg/ml), and highly transient ( $t_{1/2}$  = 5-7 minutes). To circumvent this problem, it is necessary to measure a metabolite that cannot be formed by platelets or by the kidney. TXB<sub>2</sub> can be metabolized by 11-hydroxy TX dehydrogenase to form 11-dehydro TXB<sub>2</sub>, or by β-oxidation to form 2.3-dinor TXB<sub>3</sub>. Infusion studies using TXB<sub>3</sub> have shown that both metabolites are formed equally, although 11-dehydro TXB, has a longer circulating half-life ( $t_{1/2}$  = 45 minutes). Therefore, measurement of 11-dehydro TXB<sub>2</sub> in plasma or urine will give a time-integrated indication of TXA<sub>2</sub> production.

96 strip/solid wells 480 strip/solid wells



#### 2.3-dinor Thromboxane B<sub>2</sub>

[63250-09-9]

MF:  $C_{18}H_{30}O_6$  FW: 342.4 Purity:  $\geq 98\%^*$ 

A solution in methyl acetate **Stability:** ≥1 year at -20°C

Summary: An abundant urinary metabolite of TXB2 that can be used as a marker for TXA2 synthesis in vivo

19500 2,3-dinor Thromboxane B<sub>2</sub> EIA Kit

519501

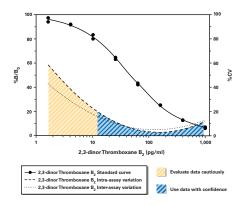
**Stability:** ≥1 year at -20°C

Sensitivity: 50% B/B $_0$ : 35 pg/ml • 80% B/B $_0$ : 7 pg/ml

Summary: TXA2 is produced from arachidonic acid by many cells and causes irreversible platelet aggregation and vascular and bronchial smooth muscle contraction. TXA<sub>2</sub> is rapidly hydrolyzed non-enzymatically to form TXB<sub>2</sub>. Although it is common to estimate TXA2 levels by measuring TXB2, most of the TXB2 measured is due to ex vivo platelet activation or intra-renal production. Measurement errors are compounded by the fact that normal concentrations of circulating TXB<sub>2</sub> are extremely low (1-2 pg/ml), and highly transient ( $t_{1/2} = 5-7$  minutes). To circumvent this problem, it is necessary to measure a metabolite that cannot be formed by platelets or by the kidney. TXB<sub>2</sub> may be metabolized by 11-hydroxy TX dehydrogenase to form 11-dehydro TXB<sub>2</sub>, or by  $\beta$ -oxidation to form 2,3-dinor TXB<sub>2</sub>. Infusion studies using TXB<sub>2</sub> have shown that both metabolites are formed equally, but that the circulating half-life of 2,3-dinor TXB<sub>2</sub> is shorter ( $t_{1/2} = 15$ minutes). Therefore, measurement of 2,3-dinor TXB2 will give a more episodic indication of TXA2 production.

519051

96 strip/solid wells 480 strip/solid wells



#### Thromboxane B<sub>2</sub> $[71953-80-5] \Delta^{17}-TXB_2$

MF:  $C_{20}H_{32}O_6$  FW: 368.5 Purity:  $\geq 98\%^*$ 

A solution in methyl acetate **Stability:** ≥2 years at -20°C

Summary: Stable hydrolysis product of TXA<sub>2</sub> synthesized from EPA by COX and TX synthase

50 µg 100 µg 500 µg 1 mg

19050

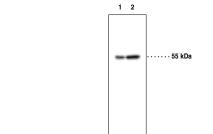
19990

• Also Available: 11-dehydro Thromboxane B<sub>3</sub> (19995)

#### Thromboxane Synthase Polyclonal Antibody 160715

Affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human thromboxane synthase amino acids 359-377 • Host: rabbit • Cross Reactivity: (+) human, porcine, murine, and rat TX synthase • Application(s): IHC and WB • TX synthase catalyzes the conversion of PGH2 to TXA2, which is a potent vasoconstrictor and inducer of platelet aggregation.



\*All 5-cis 2-series PGs (those containing a 5,6-double bond) will contain a small amount of the 5-trans ison This isomer is generally undetectable using normal phase silica columns and plates, but may be resolved using RP-HPLC. The purity for all such 2-series PGs excludes the 1-3% trans isomer which will generally be present • Also Available: Thromboxane Synthase Blocking Peptide (360715) [54857-86-2] RMI 14514, 5-(Tetradecyloxy)-2-Furoic Acid

MF:  $C_{19}H_{32}O_4$  FW: 324.5 Purity:  $\geq 98\%$ 

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: An inhibitor of fatty acid synthesis that blocks the synthesis of malonyl-CoA by acetyl-CoA carboxylase

10 mg 50 mg 100 mg

#### TP Receptor (human) Polyclonal Antibody 10004452

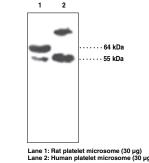
Thromboxane A2 Receptor

Peptide affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: human TP receptor C-terminal amino acids 323-343 • Host: rabbit • Cross Reactivity: (+) human, murine, rat, and Cos-7 (African green monkey) TP receptor • Application(s): ICC and WB • The TP receptor is a GPCR that mediates the action of TXA2.

500 µl

**TOFA** 



• Also Available: TP Receptor (human) Blocking Peptide (10009368)

#### TP Receptor (human) Polyclonal FITC Antibody

Thromboxane A, Receptor

Fluorescein-labeled peptide affinity-purified IgG Stability: ≥1 year at -20°C

Summary: Antigen: human TP receptor C-terminal amino acids 323-343 • Host: rabbit • Cross Reactivity: (+) Human, murine, rat, and Cos-7 (African green monkey) TP receptor • Application(s): FC, IF, and WB • The TP receptor is a GPCR that mediates the action of TXA<sub>2</sub>.

500 µl

#### TP Receptor (murine) Polyclonal Antibody 101882

Thromboxane A, Receptor

Affinity-purified IgG **Stability:** ≥1 year at -20°C

Summary: Antigen: murine TP receptor amino acids 275-289 • Host: rabbit • Cross Reactivity: (+) human, murine, rat, Cos-7 (African green monkey), and bovine TP receptor (able to detect both TPα and TPβ isoforms) • Application(s): IHC and WB • The TP receptor is a GPCR that mediates the action of TXA<sub>2</sub>.

• Also Available: TP Receptor (murine) Blocking Peptide (10004110)

#### 10162 Treprostinil

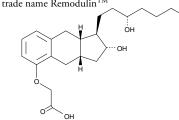
[81846-19-7]

**MF:** C<sub>23</sub>H<sub>34</sub>O<sub>5</sub> **FW:** 390.5 **Purity:** ≥98%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A stable analog of PGI, that is used clinically for the treatment of primary pulmonary hypertension under the trade name Remodulin<sup>TM</sup>

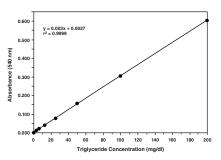
1 mg 5 mg 10 mg 50 mg



#### 10005263 Trialyceride Assay Kit

**Stability:** ≥6 months at -20°C

Summary: The measurement of TG levels are useful in the diagnosis of primary and secondary hyperlipoproteinemia, dyslipidemia, and triglyceridemia. Cayman's TG Assay provides a simple, reproducible, and sensitive tool for measuring TGs in plasma and serum. The assay is initiated with the enzymatic hydrolysis of TGs by lipase to produce glycerol and free fatty acids. The glycerol released is subsequently measured by a coupled enzymatic reaction system with a colorimetric readout at 540 nm.



#### U-18666A

10009085

[3039-71-2]

10012559

**MF:** C<sub>24</sub>H<sub>41</sub>NO<sub>2</sub> • HCl **FW:** 412.1 **Purity:** ≥95%

A crystalline solid **Stability:** ≥2 years at -20°C

Summary: A cell permeable drug that inhibits cholesterol trafficking from late endosomes/lysosomes to the ER, but not to the plasma membrane

5 mg 10 mg 25 mg 50 mg

#### U-44069 16440

[56985-32-1] 9,11-epoxymethano PGH $_{2}$ , 9,11-dideoxy-9lpha,11lpha-epoxymethano PGF $_{2lpha}$ MF:  $C_{21}H_{34}O_4$  FW: 350.5 Purity:  $\geq 98\%^*$ 

A solution in methyl acetate **Stability:** ≥2 years at -20°C

Summary: A stable analog of PGH2 and a TP receptor agonist; stimulates shape change in human platelets without a measurable increase in Ca2+ with an EC50 value of 1.8 nM

1 mg 5 mg 10 mg 50 ma

#### [56985-40-1] 9,11-dideoxy- $9\alpha$ ,11 $\alpha$ -methanoepoxy PGF<sub>2 $\alpha$ </sub>

MF:  $C_{21}H_{34}O_4$  FW: 350.5 Purity:  $\geq$ 98%\*

A solution in methyl acetate **Stability:** ≥2 years at -20°C

Summary: A stable analog of PGH2 and a TP receptor agonist; exhibits EC50 values of 4.8, 6.0, and 7.3 nM for shape change in human, rat, and rabbit platelets, respectively, and 82, 145, and 65 nM for aggregation, respectively

1 mg 5 mg 10 mg 50 mg

16450

46 Cayman Chemical caymanchem.com INDICES

Inhibitors of Lipoprotein Modifying Enzymes				
Item No.	Item Name	Target Enzyme	IC <sub>50</sub> values	Sizes
10006482	CAY10485	ACAT-1 and 2	96 μM ACAT-1; 81 μM ACAT-2	5 mg • 10 mg • 50 mg • 100 mg
10006452	CAY10486	ACAT-1 and 2	60 μM (both enzymes)	5 mg • 10 mg • 25 mg • 50 mg
10007875	CAY10499	Hormone Sensitive Lipase	90 nM (human)	1 mg • 5 mg • 10 mg • 25 mg
10006782	Oleic Acid-2,6-diisopropylanilide	ACAT	7 nM	5 mg • 10 mg • 50 mg • 100 mg
10006529	Oleyl Anilide	ACAT	26 μΜ	5 mg • 10 mg • 50 mg • 100 mg

For a full product listing, please visit www.caymanchem.com

 $\Delta^{17}$ -U-46619

1646

16465

MF:  $C_{21}H_{32}O_4$  FW: 348.5 Purity:  $\geq$ 98%\* A solution in methyl acetate **Stability:**  $\geq$ 2 years at -20°C **Summary:** A stable analog of TXA<sub>3</sub> and U-46619

50 μg 100 μg 500 μg 1 mg

U-51605
[64192-56-9]

**MF:**  $C_{20}H_{32}N_2O_2$  **FW:** 332.5 **Purity:**  $\geq$ 98%\*

A solution in methyl acetate **Stability:** ≥1 year at -20°C

**Summary:** A stable analog of  $PGH_2$  that acts as an inhibitor of both PGI and TX synthases

100 µg 500 µg 1 mg 5 mg

Vasoactive Eicosanoid HPLC Mixture

Purity: ≥98% for each compound

A solution in methyl acetate **Stability:**  $\geq$ 6 months at -20°C **Summary:** Contains TXB<sub>2</sub>, 11-dehydro TXB<sub>2</sub>, 6-keto PGF<sub>1 $\alpha$ </sub>, 2,3-dinor-6-keto PGF<sub>1 $\alpha$ </sub> (100  $\mu$ g each), and 12(S)-HHTrE (5  $\mu$ g)

1 ea

(±)-Warfarin

13566

10003

[81-81-2] Athrombine-K, Coumadin, Coumafene, WARF 42, Zoocoumarin

**MF:**  $C_{19}H_{16}O_4$  **FW:** 308.3 **Purity:**  $\geq$ 98%

A crystalline solid **Stability:** ≥2 years at -20°C

**Summary:** An anti-coagulant used to prevent heart attacks, stroke, and the formation of blood clots; also used as a rodenticide

100 mg 250 mg 500 mg 1 g

• Also Available: (+)-Warfarin (13526) (-)-Warfarin (13531)

# \*All 5-cis 2-series PGs (those containing a 5,6-double bond) will contain a small amount of the 5-trans isomer. This isomer is generally undetectable using normal phase silica columns and plates, but may be resolved using RP-HPLC. The purity for all such 2-series PGs excludes the 1-3% trans isomer which will generally be present.

# Alphabetical Index

4-922500	
ACAT-1 Blocking Peptide	
ACAT-2 Blocking Peptide	
ACAT-1 Polyclonal Antibody	
ACAT-2 Polyclonal AntibodyACENOCOUNTIES (ACENOCOUNTIES )	6
Acenocoumarin (Acenocoumaroi)	
R)-(+)-Acenocoumarol	
S)-(-)-Acenocoumarol	6
Acetyl Lysine Monoclonal Antibody (Clone 7F8)	27
Acetyl Podocarpic Acid Anhydride	6
Acetylsalicyclic Acid Matol ester (Aspalatone)	8
Acetylsalicyclic Acid (Aspirin)	8
N-Acetyl Ser-Asp-Lys-Pro (AcSDKP EIA Kit)	6
ACSDKP EIA Kit	6
ACT-293987 (NS-304) Acyl-coenzyme A:Cholesterol Acyltransferase-1 (ACAT-1)	30
Acyl-coenzyme A:Cholesterol Acyltransferase-1 (ACAT-1) Acyl-coenzyme A:Cholesterol Acyltransferase-2 (ACAT-2)	
N-(5-adamantane-1-yl-methoxy-pentyl)-Deoxynojirimycin)	0
AMP-Deoxynojirimycin)	7
Adamantane-pentyl-dNM (AMP-Deoxynojirimycin)	7
ADH (Arginine Vasopressin EIA Kit)	8
Adipose Triglyceride Blocking Peptide	7
Adipose Triglyceride Lipase Polyclonal Antibody	
AM3102	34
AMP-Deoxynojirimycin	7
AMP-dNM (AMP-Deoxynojirimycin)	
Angiotensin II EIA KitAnthropodeoxycholic Acid)	/
Antidiuretic Hormone (Arginine Vasopressin EIA Kit)	۱4
APD (Acetyl Podocarpic Acid Anhydride)	٥
Apelin-13	7
Annal Blocking Pentide	7
ApoAl Monoclonal Antibody (Clone CC3821C4)	7
ApoAl Polyclonal Antibody	7
ApoAl Western Ready Control	
Apolipoprotein Al	See ApoA <u>l</u>
ATGL (Adipose Triglyceride Lipase Polyclonal)	7
20-carboxy Arachidonic Acid	
o-3 Arachidonic Acid Arginine Vasopressin EIA Kit	
Argipressin (Arginine Vasopressin EIA Kit)	0
	×
Argipressin (Arginine vasopressin EIA NI)	88 8
Aspalatone	8
Aspalatone	8 8
Aspalatone	8 8 8
Aspalatone	8 8 8 8
Aspalatone	
Aspalatone	
Aspalatone Aspirin Aspirin MEffect-Detection Kit	
Aspalatone	
Aspirin   Aspirin   Aspirin   Aspirin   Effect-Detection Kit   Atherosclerosis Product Pack   Athropeptin (rat) EIA Kit   Athrombine-K ((±)-Warfarin)   AUDA   AVP (Arginine Vasopressin EIA Kit)   Azelaoyl PAF   BAY-41-8543   BBR (Berberine)   Benzofibrate (Bezafibrate)   I-Benzylimidazole   Beraprost (sodium salt)   Berberine   Bezafibrate   Bezafibrate   Bezafibrate   Bezafibrate   Bezafibr (Bezafibrate)   Bezaflor (Bezafibrate)   BESESTERON   BESE	
Aspalatone	
Aspalatone Aspirin Aspirin Meffect-Detection Kit	
Aspalatone  Aspirin  Aspirin  Afherosclerosis Product Pack  Atriopeptin (rat) EIA Kit  Athrombine-K ((±)-Warfarin)  AUDA  AVP (Arginine Vasopressin EIA Kit)  Azelaoyl PAF  BAY-41-8543  BBR (Berberine)  Benzofibrate (Bezafibrate)  Beraprost (sodium salt)  Berberine  Bezafibrate  Bezafibrate  Bezafibrate  Bezafibrate  Bezafibrate  Bezafibrate  Bezafibrate  Bezafibrote  Botoprofic Bezafibrate)  Butanoyl PAF	
Aspalatone  Aspirin  Aspirin  Aspirin  Aspirin  Aspirin  Aspirin  Effect-Detection Kit	
Aspalatone Aspirin Aspirin Meffect-Detection Kit	
Aspalatone	
Aspalatone Aspirin Aspirin Aspirin Afherosclerosis Product Pack Atriopeptin (rat) EIA Kit Athrombine-K ((±)-Warfarin) AUDA AVP (Arginine Vasopressin EIA Kit) Azelaoyl PAF BAY-41-8543 BBR (Berberine) Benzofibrate (Bezafibrate) Benzofibrate (Bezafibrate) Besports (sodium salt) Bescalip (Bezafibrate) Bothology (Bezafib	
Aspirin Maspirin Effect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspirin Maspirin Effect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspalatone Aspirin Aspirin Aspirin Aftherosclerosis Product Pack Atriopeptin (rat) EIA Kit Athrombine-K ((±)-Warfarin) AUDA AVP (Arginine Vasopressin EIA Kit) Azelaoyl PAF Asy-41-8543 BBR (Berberine) Benzofibrate (Bezafibrate) Benzofibrate (Bezafibrate) Berberine Bezafibrate Bezafi	
Aspirin Meffect-Detection Kit	
Aspirin Maspirin Effect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspirin Meffect-Detection Kit	
Aspirin Meffect-Detection Kit	

CAY1049911,46	
CAY1051411,34	
CAY1053521,23	
CAY1057334	
CAY10591	
CAY10592	
CAY10603 27	
CBHA	
CD36 Blocking Peptide	
CD36 Monoclonal Antibody (Clone JC63.1)11	
CD36 Monoclonal Antibody (Clone JC63.1) (azide free)11	
CD36 Monoclonal FITC Antibody (Clone JC63.1)11	l
CD36 Polyclonal Antibody	1
CDCA (Chenodeoxycholic Acid)	ł
Cetaben	
CG 4203 (Taprostene (sodium salt))	
CG 4305	
Chenodeoxycholic Acid14	1
Chidamide	7
Cholestane-6-oxo-3 $\beta$ ,5 $\alpha$ -diol (5 $\alpha$ -hydroxy-6-keto Cholesterol)15	5
Cholesterol Assay Kit	
Cholesterol Cell-Based Detection Assay Kit	
3-dodecanoyl-NBD Cholesterol	
3-hexanoyl-NBD Cholesterol	
Cholesteryl Linoleate Hydroperoxides	Ś
Cicaprost	5
Ciglitazone	
Ciloprost (Iloprost)	2
Ciprostene (calcium salt)	,
15,34	
(±)-Clopidogrel (hydrochloride)	)
3-hydroxy-3-methylglutaryl-Coenzyme A-d <sub>3</sub> (ammonium salt)	)
20-COOH-AA (20-carboxy Arachidonic Acid)	7
Coumadin ((±)-Warfarin)	5
Coumafene ((±)-Warfarin)	Ś
coumarin-SAHA27	7
COX-1 Monoclonal Antibody (Clone CX111)16	
COX-1 Monoclonal FITC Antibody (Clone CX111)	5
COX-1 Monoclonal PE Antibody (Clone CX111)	)
COX-1 (minite) Polyclonal Antibody	
COX-2 Monoclonal Antibody (Clone CX229)16	5
COX-2 Monoclonal FITC Antibody (Clone CX229)	5
COX-2 (human) Blocking Peptide	
COX-2 (human) Polyclonal Antibody	ò
Goat Anti-COX-2 (human) Affinity-Purified Polyclonal Antibody	(
COX-2 (murine) Polyclonal Antibody	
COX-2 (murine) Polyclonal FITC Antibody	Ś
COX-2 (murine) Polyclonal Antiserum	Ś
cPGI (Carbaprostacyclin)10	
C-Reactive Protein (human) EIA Kit10	)
CRP (C-Reactive Protein (human) EIA Kit))	
CS 500 (Mevastatin)	5
CUDA	
Cyclooxygenase 1 (COX-1 Monoclonal Antibody (Clone CX111))	
Cyclooxygenase 2 (COX-2 Monoclonal Antibody (Clone CX229))	
(1R,2R)-2-(4'-(3-phenylureido)biphenylcarbonyl)-cyclopentanecarboxylic	
acid) (A-922500)6	
Cytosolic Epoxide Hydrolase See Soluble Epoxide Hydrolase	
DCU (N,N'-Dicyclohexylurea)	
11-dehydro Thromboxane B <sub>2</sub> 44	1
11-dehydro Thromboxane B <sub>2</sub> d <sub>4</sub>	1
11-dehydro Thromboxane B <sub>2</sub> EIA Kit44	4
11-dehydro Thromboxane Ballipid Maps MS Standard 44	1
11-dehydro Thromboxane B <sub>2</sub> Quant-PAK	1
11-dehvdro Thromboxane B <sub>2</sub> 44	ļ
N-(5-adamantane-1-yl-methoxy-pentyl)-Deoxynojirimycin) (AMP-	7
Deoxynojirimycin)	1
Desnutrin (Adipose Triglyceride Lipase Polyclonal Antibody)	7
N N'-Dicyclohexylured	4
9,11-dideoxy- $9\alpha$ ,11 $\alpha$ -epoxymethano PGF <sub>2<math>\alpha</math></sub> (U-44069)	5
9,11-dideoxy-9α,11α-methanoepoxy PGF <sub>2a</sub> (U-44619)	5
Difaterol (Bezafibrate)	
5(\$),6(R)-DiHETE	
5(S),6(R)-DIHETE Lipid Maps MS Standard	
5(S) -6(S) -DIHETE	1
5(S),6(S)-DiHETE Lipid Maps MS Standard41	ı
5(\$),15(\$)-DiHETE41	ı

E(C) 1 E(C) D'HETE L'. '.   1 4 1 (C)
5(S),15(S)-DiHETE Lipid Maps MS Standard41
8(S),15(S)-DiHETE
(±)14,15-DiHETE
3,4-dihydroxy Hydrocinnamic Acid (L-Aspartic Acid dibenzyl ester) amide
(CAY10485)
3,4-Dihydrocinnamic Acid (L-alanine methyl ester) amide (CAY10487)11
2.3-dinor-6-keto Prostaglandin F. FIA Kit 38
2,3-dinor-6-keto Prostaglandin $F_{1\alpha}$ (sodium salt)
2,3-dinor-6-keto Prostaglandin $F_{1\alpha}$ -d <sub>4</sub> (sodium salt)
2,3-dinor Thromboxane B <sub>2</sub> EIA Kit
Diphenyl-1-pyrenylphosphine
Docosahexaenoic Acid
Docosapentaenoic Acid
Docosatrienoic Acid
3-dodecanoyl-NBD Cholesterol14
DPPP (Diphenyl-1-Pyrenylphosphine)
DuP 753 (Losartan (potassium salt))
ecNOS
14,15-EE-5(Z)-E
14,15-EE-8(Z)-E
(±)5(6)-EET40
(±)5(6)-EET-d <sub>11</sub> 40
(±)5(6)-EET Ethanolamide40
(±)5(6)-EET Lipid Maps MS Standard40
(±)5(6)-EET methyl ester
(±)8(9)-EET
(±)8(9)-EE1-a <sub>11</sub>
(±)8(9)-EET Lipid Maps MS Standard
(±)8(9)-EET methyl ester
(±)11(12)-EET
(±)11(12)-EET-d <sub>11</sub>
(±)11(12)-FET Ethanolamide40
(±)11(12)-EET Lipid Maps MS Standard40
(±)11(12)-EET methyl ester40
(±)14(15)-EET40
(±)14(15)-EET-d <sub>11</sub> 40
(±)14(15)-EET Ethanolamide
(±)14(15)-EET Lipid Maps MS Standard
(±)14(15)-EET methyl ester
(±) 14(13)-EE1-31
Eicosapentaenoic Acid-d <sub>s</sub>
Eicosapentaenoic Acid ethyl ester
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelia ElA Kit 17 eNOS Blocking Peptide 29 eNOS (bovine recombinant) 29 eNOS (Electrophoresis Standard 29 eNOS Polyclonal Antiserum 29
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (Electrophoresis Standard 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 4(5)-EpDPE methyl ester 40 19(20)-EpDPE 40
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelia ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS Electrophoresis Standard 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 4(5)-EpDPE methyl ester 40 16(17)-EpDPE 40 19(20)-EpDPE 40 5(6)-EpETE methyl ester 40
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free)       17         EL (Endothelial Lipase Polyclonal Antibody)       17         Endothelial (human) Lipase Blocking Peptide       17         Endothelial Lipase Polyclonal Antibody       17         Endothelia ElA Kit       17         eNOS Blocking Peptide       29         eNOS (bovine recombinant)       29         eNOS (bectrophoresis Standard       29         eNOS Polyclonal Antiserum       29         EPA-EE (Eicosapentaenoic Acid ethyl ester)       EPA ethyl ester         (Eicosapentaenoic Acid ethyl ester)       17         4(5)-EpDPE methyl ester       40         19(20)-EpDPE       40         19(20)-EpDPE       40         15(4)5-EpETE methyl ester       40         14(15)-EpETE       40         17(18)-EpETE       40
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (bovine recombinant) 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 ENDPE methyl ester 40 EPA-EDPE Methyl ester 40 EPA-EE (Eicosapentaenoic Acid ethyl ester) 40 EPA-EE (Eicosapentaenoic Acid
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (Electrophoresis Standard 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 4(5)-EpDPE methyl ester 40 19(20)-EpDPE 40 19(20)-EpDPE 40 19(15)-EpETE 40 14(15)-EpETE 40 14(15)-EpETE 40 14(15)-EpETE 40 14(15)-EpETE 40 14(15)-EpETE 40 15(6)-EpETE 50 50 Juble Epoxide Hydrolase (±)-9(10)-EpOME 40
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelia ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS Electrophoresis Standard 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 EATO EPA ETHYL EPA ETHYL ENDE 17 ENDE methyl ester 17 ENDE methyl ester 17 ENDE 17
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (bovine recombinant) 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 ENGS Electrophoresis Standard 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 17 EPA ethyl ester 17 EPA ethyl ester 17 EPA ethyl ester 19 EPA ethyl ester 19 EPA ethyl ester 19 EPA ethyl ester 19 EPA ethyl ester 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 40 EP
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (bovine recombinant) 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 ENGS Electrophoresis Standard 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 17 EPA ethyl ester 17 EPA ethyl ester 17 EPA ethyl ester 19 EPA ethyl ester 19 EPA ethyl ester 19 EPA ethyl ester 19 EPA ethyl ester 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 40 EP
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (bovine recombinant) 29 ENOS Electrophoresis Standard 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 17 Elicosapentaenoic Acid ethyl ester) 17 EPA ethyl ester 17 EPA ethyl ester 17 EPA ethyl ester 17 EPA ethyl ester 19 EPA ethyl es
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (bovine recombinant) 29 ENOS Electrophoresis Standard 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 17 Elicosapentaenoic Acid ethyl ester) 17 Elicosapentaenoic Acid ethyl ester) 17 EPA ethyl ester 17 Elicosapentaenoic Acid ethyl ester) 17 EPA ethyl ester 17 EPA ethyl
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelia ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS Electrophoresis Standard 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 ENDET methyl ester 17 ENDET methyl ester 18 EPA ethyl ester 19 EPA
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENO'S Blocking Peptide 29 ENO'S (bovine recombinant) 29 ENO'S (bovine recombinant) 29 ENO'S Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 END'E methyl ester 17 END'E methyl ester 19 EPA-EDPE 19 EPA-
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 Endothelial ElA Kit 17 Endothelial ElA Kit 17 Endos Blocking Peptide 17 Endothelial ElA Kit 17 Endos Blocking Peptide 19 ENOS (bovine recombinant) 19 ENOS Electrophoresis Standard 19 ENOS Polyclonal Antiserum 19 EPA-EE (Eicosapentaenoic Acid ethyl ester) 19 EPA-EE (Eicosapentaenoic Acid ethyl ester) 19 EPA-EE (Eicosapentaenoic Acid ethyl ester) 19 EPA-EDPE methyl ester 19 EPA-EDPE 19 EPA-EDPE 19 EPA-EDPE 19 EPA-EDPE 19 EPA-EDPE 19 EPA-ED ElE Methyl ester 19 EPA-EB ElE
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS (bovine recombinant) 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) 17 Elosapentaenoic Acid ethyl ester) 17 Elosapentaenoic Acid ethyl ester) 17 Elosapentaenoic Acid ethyl ester) 17 EPA ethyl ester 17 EPA ethyl ester 18 EPA ethyl ester 19 EPA
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free)
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 (4(5)-EpDPE methyl ester 40 16(17)-EpDPE 40 16(17)-EpDPE 40 16(17)-EpDPE 40 16(17)-EpDPE 40 17(18)-EpETE methyl ester 40 17(18)-EpETE 40 17(18)-EpETE 40 17(18)-EpETE 40 17(18)-EpETE 40 17(18)-EpOME 40 (±)-9(10)-EpOME-44 40 (±)-12(13)-EpOME-44 40 (±)-1
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial Lipase Polyclonal Antibody 17 Endothelial ElA Kit 17 Endothelia Ela Kit 18 Ela Kit 19 Endothelia Ela Kit 19 Endothe
Eicosapentaenoic Acid (peroxide free) 17 EL (Endothelial Lipase Polyclonal Antibody) 17 Endothelial (human) Lipase Blocking Peptide 17 Endothelial Lipase Polyclonal Antibody 17 ENOS Blocking Peptide 29 ENOS (bovine recombinant) 29 ENOS Polyclonal Antiserum 29 EPA-EE (Eicosapentaenoic Acid ethyl ester) EPA ethyl ester (Eicosapentaenoic Acid ethyl ester) 17 (4(5)-EpDPE methyl ester 40 16(17)-EpDPE 40 16(17)-EpDPE 40 16(17)-EpDPE 40 16(17)-EpDPE 40 17(18)-EpETE methyl ester 40 17(18)-EpETE 40 17(18)-EpETE 40 17(18)-EpETE 40 17(18)-EpETE 40 17(18)-EpOME 40 (±)-9(10)-EpOME-44 40 (±)-12(13)-EpOME-44 40 (±)-1

5-(Tetradecyloxy)-2-Furoic Acid (TOFA)	45
G-23350 (Acenocoumarol) Goat Anti-COX-2 (human) Affinity-Purified Polyclonal Antibody	6
GPIIIbGPIIIb	16
GPIV	
GSK264220A	18
Guanylate Cyclase $lpha_1$ subunit (soluble) Blocking Peptide Guanylate Cyclase $lpha$ subunit (soluble) Polyclonal Antibody	
Guanylate Cyclase β1 subunit (soluble) Blocking Peptide	18
Guanylate Cyclase β, subunit (soluble) Polyclonal Antibody	18
E)-Guggulsterone	18
2)-Guggusterone	18.34
GW 7647	19,34
GW 9578	
GW 9662	
HAT Inhibitor Screening Assay Kit	
HDAC Activity Assay Kit	
HDAC Cell-Based Activity Assay Kit	2/
HDAC1 Inhibitor Screening Assay Kit	27
HDAC1 Polyclonal Antibody	27
HDAC2 (human recombinant) HDAC3/NCOR2 (human recombinant)	27
HDAC3 Polyclonal Antibody	
HDAC4 (human recombinant)	27
HDAC4 Polyclonal Antibody	
HDAC5 (human recombinant)HDAC6 (human recombinant)	2/
HDAC6 Polyclonal Antibody	27
HDAC7 (Phospho-Ser <sup>155</sup> ) Polyclonal Antibody	27
HDAC8 (human recombinant)	27
HDAC9 (human recombinant)	27
HDAC11 Polyclonal Antibody	27
S)-HDAC-42	27
Hekbilin (Chenodeoxycholic Acid)	14
HET0016	
19(R)-HETE	
19(S)-HETE P-cis-Hexadecenoic Acid (Palmitoleic Acid)	
Hexadecvlamino-p-amino Benzoic Acid (Cetaben)	11
3-hexanoyl-NBD Cholesterol	14
Hexarelin ReceptorHistone H2B (Xenopus recombinant)	
Historie HZB (Neriopus recombinant; G103A subsitution)	27
Histone H4 (human recombinant)	27
HMG-CoA-d <sub>3</sub> (3-hydroxy-3-methylglutaryl- Coenzyme A-d <sub>3</sub> (ammonium salt))	1.5
HNHA	27
±)9-HODE cholesteryl ester	19
P(R)-HODE cholesteryl ester	
P(S)-HODE cholesteryl ester±)13-HODE cholesteryl ester	
13(R)-HODE cholesteryl ester	
13(S)-HODE cholesteryl ester	19
Hormone Sensitive Lipase (human recombinant)	22
Hormone Sensitive Lipase Polyclonal Antibody	
HSLSee Hormone Sensi	
HTS 01037	
3,4-dihydroxy Hydrocinnamic Acid (L-Aspartic Acid dibenzyl ester) (CAY10485)	amide
N-hydroxy-N'-(4-n-butyl-2-methylphenyl)Formamidine (HET0016)	19
1-Hydroxycinnamic Acid (L-phenylalanine methyl ester)	
amide (CAY10486)	11
5α-hydroxy-6-keto Cholesterol 3-hydroxy-3-methylglutaryl-Coenzyme A-d <sub>3</sub> (ammonium salt)	
-BOP	22,23
loprost	
odophenyl sulfonyl amino pinane TXA <sub>2</sub> (I-SAP)	
4-iodo-SAHAP Receptor (human) Blocking Peptide	
P Receptor (human) Polyclonal Antibody	22
P Receptor (murine) Blocking Peptide	22
P Receptor (murine) Polyclonal AntibodySAP	
KB2115	23
(DdiA-PC	23
Kebilis (Chenodeoxycholic Acid)	
S-keto Prostaglandin F $_{1lpha}$ S-keto Prostaglandin F $_{1lpha}$ -d $_4$	აგ 38
6-keto Prostaglandin F <sub>12</sub> -d <sub>4</sub> Lipid Maps MS Standard	38
S-keto Prostaglandin F <sub>1α</sub> EIA Kit	38

	3
11-keto TXB <sub>2</sub> (11-dehydro Thromboxane B <sub>2</sub> )	4
KOdiA-PC	2
- 637312 (Mevastatin)	2
655,240	
CAT Blocking Peptide	2
CAT Polyclonal Antibody	2
DL Receptor Blocking Peptide	
DL Receptor Polyclonal Antibody	2
DLR (LDL Recptor Polyclonal Antibody)	
DL Uptake Cell-Based Assay Kit	2
ecithin:Cholesterol Acyltransferase (LCAT Polyclonal Antib	oody)2
eukotriene A <sub>4</sub> Hydrolase (human recombinant)	2
inolein Hydroperoxides	2
x-Linolenić Acid	1
ipid Droplets Fluorescence Assay Kit	2
ipoprotein PLA <sub>2</sub> See PA	AF Acetylhydrolase
5-Lipoxygenase Blocking Peptide	2
5-Lipoxygenase (human recombinant)	
5-Lipoxygenase Polyclonal Antibody	2
5-Lipoxygenase (Phospho-Ser523) Polyclonal Antibody	2
2-Lipoxygenase (murine leukocyte) Polyclonal Antiserum	2
2-Lipoxygenase (platelet-type, murine recombinant)	2
5-Lipoxygenase-1 (rabbit) Polyclonal Antiserum	2
isinopril	
osartan (potassium salt)	2
ovastatin	14,2
.ovastatin Hydroxy Acid (sodium salt)	14,2
p-PLA <sub>2</sub> See PA	AF Acetylhydrolasi
TA,H (Leukotriene A, Hydrolase (human recombinant))	2
uminex® Thromboxane B. Kit	2
uminex <sup>®</sup> 11-dehydro Thromboxane B <sub>2</sub> Kit	2
yso-PAF C-16	3
yso-PAF C-16-d <sub>4</sub>	ર
yso-PAF C-18	
yso-PAF C-18-d <sub>4</sub>	
pxLDL Receptor	
A 344	
MEDCIA 16	
0.11 mathana anovy Prostaglandin E	رح
$9,11$ -methane-epoxy Prostaglandin $F_{1\alpha}$ $9,11$ -dideoxy- $9\alpha,11\alpha$ -methanoepoxy PG $F_{2\alpha}$ (U-44619)	د
/othyl 2 bydrovy 2 /2 poptyl ovyphonyl) oct 5 ypogto (C/	4. V10514\ 1
Methyl-8-hydroxy-8-(2-pentyl-oxyphenyl)-oct-5-ynoate (CA	1110514)1
2-O-methyl PAF C-16	د
2-O-methyl PAF C-18	
Mevastatin	
AL 236B (Mevastatin)	
MK 521 (Lisinopril)	2
MK 522 (Lisinopril)	2
MK 733 (Simvastatin)	4
MK 954 (Losartan (potassium salt))	2
AL 1129 (Beraprost (sodium salt))	
MRE-269	
AS-275	
Myeloperoxidase Chlorination Assay Kit	
Myeloperoxidase (human) EIA Kit	2
Myeloperoxidase Inhibitor Screening Assay Kit	2
Myeloperoxidase Peroxidation Assay Kit	2
Ayricetin	2
VÁRC-1	See PCSK
-NEPC	2
licoumalone (Acenocoumarol)	2
licoumalone (Acenocoumarol) IOS III	
licoumalone (Acenocoumarol)	
licoumalone (Acenocoumarol)	
licoumalone (Acenocoumarol) IOS III IS-304	
licoumalone (Acenocoumarol) IOS III IS-304	
licoumalone (Acenocoumarol) IOS III IS-304 ISC 17013 (N,N'-Dicyclohexylurea) ISC 281245 (Mevastatin) ISC 354317 (Resorcinonaphthalein)	
licoumalone (Acenocoumarol)	
licoumalone (Acenocoumarol)	
licoumalone (Acenocoumarol) IOS III IS-304 ISC 17013 (N,N'-Dicyclohexylurea). ISC 281245 (Mevastatin). ISC 354317 (Resorcinonaphthalein). ISC 407290 (Myricetin)	
licoumalone (Acenocoumarol) IOS III IS-304 ISC 17013 (N,N'-Dicyclohexylurea) ISC 281245 (Mevastatin) ISC 354317 (Resorcinonaphthalein) ISC 407290 (Myricetin) IUC 1 Iuclear Hormone Receptor 1 Iucley Anilide) I-Octadecyl-N'-propyl sulfamide	
licoumalone (Acenocoumarol) IOS III IS-304 ISC 17013 (N,N'-Dicyclohexylurea) ISC 281245 (Mevastatin) ISC 354317 (Resorcinonaphthalein) ISC 407290 (Myricetin) IUC 1 Iuclear Hormone Receptor 1 Iucley Anilide) I-Octadecyl-N'-propyl sulfamide	
Aicoumalone (Acenocoumarol)  AIS-304  AIS-304  AIS-304  AISC 17013 (N,N'-Dicyclohexylurea)  AISC 281245 (Mevastatin)  AISC 354317 (Resorcinonaphthalein)  AISC 407290 (Myricetin)  AIUC 1  Auclear Hormone Receptor 1  DA (Oleyl Anilide)  AISC AISC AISC AISC AISC AISC AISC AISC	
Jicoumalone (Acenocoumarol)  JOS III  JIS-304  JISC 17013 (N,N'-Dicyclohexylurea)	
Alicoumalone (Acenocoumarol) AOS III AIS-304 AISC 17013 (N.N'-Dicyclohexylurea) AISC 281245 (Mevastatin) AISC 354317 (Resorcinonaphthalein) AISC 407290 (Myricetin) AISC 40729	
Aicoumalone (Acenocoumarol) AICS III AIS-304 AISC 17013 (N.N'-Dicyclohexylurea) AISC 281245 (Mevastatin) AISC 354317 (Resorcinonaphthalein) AISC 407290 (Myricetin) AIUC1 AIUC	
Aicoumalone (Acenocoumarol)  AIS-304  AIS-304  AISC 17013 (N,N'-Dicyclohexylurea)  AISC 281245 (Mevastatin)  AISC 354317 (Resorcinonaphthalein)  AISC 407290 (Myricetin)  A	
Nicoumalone (Acenocoumarol)  NOS III  NS-304  NSC 17013 (N,N'-Dicyclohexylurea)  NSC 281245 (Mevastatin)  NSC 354317 (Resorcinonaphthalein)  NSC 407290 (Myricetin)  NUC1  NUC1  NUClear Hormone Receptor 1  DA (Oleyl Anilide)  N-Octadecyl-N'-propyl sulfamide  DEA (Oleoyl Ethanolamide)  DKY-046 (Ozagrel)  DIEIC Acid Anilide (Oleyl Anilide)  DIEIC Acid Acid-2,6-diisopropylanilide  DIEIC Acid Ethanolamide (Oleoyl Ethanolamide)  DIEIC Acid Ethanolamide (Oleoyl Ethanolamide)	
Nicoumalone (Acenocoumarol) NOS III NOS III NSC 17013 (N,N'-Dicyclohexylurea) NSC 281245 (Mevastatin) NSC 281245 (Mevastatin) NSC 354317 (Resorcinonaphthalein) NSC 407290 (Myricetin) NUC 1 Nuclear Hormone Receptor 1 DA (Oleyl Anilide) N-Octadecyl-N'-propyl sulfamide DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) DIeic Acid Anilide (Oleyl Anilide) Dleic Acid Ethanolamide (Oleoyl Ethanolamide) Dleoyl Ethanolamide (Oleoyl Ethanolamide) Dleoyl Ethanolamide	
Nicoumalone (Acenocoumarol) NOS III NS-304 NSC 17013 (N.N'-Dicyclohexylurea) NSC 281245 (Mevastatin) NSC 354317 (Resorcinonaphthalein) NSC 407290 (Myricetin) NUC1 Nuclear Hormone Receptor 1 DA (Oleyl Anilide) N-Octadecyl-N'-propyl sulfamide DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) DEA (Odeyl Anilide (Oleyl Anilide) Dieic Acid Anilide (Oleyl Anilide) Dieic Acid Ethanolamide (Oleoyl Ethanolamide) Dieyl Anilide Dieyl Anilide Dieyl Anilide Dieyl Anilide Dixidized Lipid HPLC Mixture	
Nicoumalone (Acenocoumarol) NOS III NS-304 NSC 17013 (N.N'-Dicyclohexylurea) NSC 281245 (Mevastatin) NSC 354317 (Resorcinonaphthalein) NSC 407290 (Myricetin) NUC1 Nuclear Hormone Receptor 1 DA (Oleyl Anilide) N-Octadecyl-N'-propyl sulfamide DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) DEA (Odeyl Anilide (Oleyl Anilide) Dieic Acid Anilide (Oleyl Anilide) Dieic Acid Ethanolamide (Oleoyl Ethanolamide) Dieoyl Ethanolamide	
in NEPC.  Ilicoumalone (Acenocoumarol)  Ilicoumalone (Acenocoumarol)  Ilicoumalone (Acenocoumarol)  Ilicoumalone (Acenocoumarol)  Ilicoumalone (Acenocoumarol)  Ilicoumalone (Ilicoumarol)  Ilicoumarol	See eNO 30. 31 22 32 32 34 30,3 30,3 30,4 30,4 30,4 30,1 30,1 30,1 30,2 30,1 30,2 30,2 30,3 30,4 30,4 30,4 30,4 30,4 30,4 30,4
Nicoumalone (Acenocoumarol) NOS III NOS III NISC 17013 (N,N'-Dicyclohexylurea) NISC 281245 (Mevastatin) NISC 354317 (Resorcinonaphthalein) NISC 407290 (Myricetin) NUC1 Nuclear Hormone Receptor 1 DA (Oleyl Anilide) N-Octadecyl-N'-propyl sulfamide DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) DICIA Acid Ethanolamide (Oleoyl Ethanolamide) DICIA Acid Ethanolamide DICIA Coleoyl Ethanolamid	See eNO 3 3 3 3 See PPAR 3 3 3 3 3 3 3 3 3 3 3 3 3 3 3 3 3 3 3
Nicoumalone (Acenocoumarol) NOS III NOS III NS-304 NSC 17013 (N,N'-Dicyclohexylurea) NSC 281245 (Mevastatin) NSC 354317 (Resorcinonaphthalein) NSC 407290 (Myricetin) NUC1 NUClar Hormone Receptor 1 DA (Oleyl Anilide) DA (Oleyl Anilide) N-Octadecyl-N'-propyl sulfamide DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) DIeic Acid Anilide (Oleyl Anilide) DIeic Acid Anilide (Oleyl Anilide) DIeic Acid Ethanolamide (Oleoyl Ethanolamide) DIeyl Anilide Dieyl Anilide Dieyl Anilide Dixidized Lipid HPLC Mixture Dixidized Low-Density Lipoprotein-β₂Glycoprotein I (humarn (human) ELSA Kit) So-Oxo-3,5-diol (5α-hydroxy-6-keto Cholesterol)	
Nicoumalone (Acenocoumarol) NOS III NS-304 NSC 17013 (N.N'-Dicyclohexylurea) NSC 281245 (Mevastatin) NSC 354317 (Resorcinonaphthalein) NSC 407290 (Myricetin) NUC1 Nuclear Hormone Receptor 1 DA (Oleyl Anilide) N-Octadecyl-N'-propyl sulfamide DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) Dieic Acid Anilide (Oleyl Anilide) Dieic Acid Anilide (Oleyl Anilide) Dieic Acid Ethanolamide (Oleoyl Ethanolamide) Dieyl Anilide Dieyl Anilide Dieyl Ethanolamide Dieyl Ethanolamide Dieyl Ethanolamide Dieyl Ethanolamide Dieyl Anilide Dieyl Anilide Dieyl Sylvanion Dieyl Anilide	See eNO 3. 3. 3. 2. See PPAR See PPAR See PPAR 3. 3. 3. 3. 3. 3. 3. 3. 3. 3. 3. 3. 3.
Nicoumalone (Acenocoumarol) NOS III NOS III NS-304 NSC 17013 (N,N'-Dicyclohexylurea) NSC 281245 (Mevastatin) NSC 354317 (Resorcinonaphthalein) NSC 407290 (Myricetin) NUC1 NUClar Hormone Receptor 1 DA (Oleyl Anilide) DA (Oleyl Anilide) N-Octadecyl-N'-propyl sulfamide DEA (Oleoyl Ethanolamide) DEA (Oleoyl Ethanolamide) DIeic Acid Anilide (Oleyl Anilide) DIeic Acid Anilide (Oleyl Anilide) DIeic Acid Ethanolamide (Oleoyl Ethanolamide) DIeyl Anilide Dieyl Anilide Dieyl Anilide Dixidized Lipid HPLC Mixture Dixidized Low-Density Lipoprotein-β₂Glycoprotein I (humarn (human) ELSA Kit) So-Oxo-3,5-diol (5α-hydroxy-6-keto Cholesterol)	See eNO 30 31 22 32 See PPAR See PPAR 30,3 30,3 30,4 30,4 30,4 30,1 30,4 30,4 30,5 30,4 30,5 30,4 30,4 30,5 30,4 30,4 30,5 30,4 30,5 30,4 30,6 30,6 30,6 30,7 30,7 30,8 30,8 30,8 30,8 30,8 30,8 30,8 30,8

DAE A getulo (dralage (la unaga) Blacking Dontida	,
PAF Acetylhydrolase (human) Blocking Peptide	
PAF Acetylhydrolase (human recombinant)	
PAF Acetylhydrolase (human) Polyclonal Antibody	
PAF Acetylhydrolase (human) Western Ready Control	
PAF Acetylhydrolase Inhibitor Screening Assay Kit	
PAF-AHSee PAF	
PAF-C16	
PAF-C16-d <sub>4</sub>	
2-O-ethyl-PAF C-16	
2-O-methyl PAF C-16	
PAF C-18	
PAF C-18-d <sub>4</sub>	
2-O-methyl PAF C-18	
2-O-memyrar C-10	٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠٠
PAF Receptor Blocking Peptide (Monoclonal)	
PAF Receptor Blocking Peptide (Polyclonal)	
PAF Receptor (human) Monoclonal Antibody (11A4, Clone	21)
PAF Receptor (human) Polyclonal Antibody	
2-thio PAF	
Palmitoleate (Palmitoleic Acid)	,
Palmitoleate (Palmitoleic Acid)n-7 Palmitoleate (Palmitoleic Acid)	
Palmitalaia A aid	
Palmitoleic Acid	
cis-Palmitoleic Acid (Palmitoleic Acid)	
Pargyline (hydrochloride)	
PC (Phosphatidylcholine Assay Kit)	
pCAF Histone Acetyltransferase	
PCSK9 (human) Blocking Peptide	
PCSK9 (murine) Blocking Peptide	(
PCSK9 Monoclonal Antibody (Clone 15A6)	
PCSK9 Polyclonal Antibody	
PCSK9 (human) Polyclonal Antibody	
PCSK9 (murine) Polyclonal Antibody	
PCSK9 Western Ready Control	
Peroxisome Proliferator-Activator Receptor	See PPA
PGC-1 Blocking Peptide	33,3
PGC-1 Polyclonal Antibody	
9,11-dideoxy-9a,11a-epoxymethano PGF <sub>2a</sub> (U-44069)	
9,11-dideoxy-9a,11a-methanoepoxy PGF <sub>2a</sub> (U-44619)	
9,11-epoxymethano PGH <sub>1</sub> (9,11 methane-epoxy Prostaglan	
9,11-epoxymetriano PGH <sub>1</sub> (9,11 metriane-epoxy Prostagian	$\alpha_{l\alpha}$ /
9,11-epoxymethano PGH <sub>2</sub> (U-44069)	4
PGHS-1 (COX-1 Monoclonal Antibody (Clone CX111))	
PGHS-2 (COX-2 Polyclonal Antibody (Clone CX229))	
PGI SynthaseSee Prostac	ılandin I Synthas
PGI, Receptor	See IP Recept
PGPC.	
(1R,2R)-2-(4'-(3-phenylureido)biphenylcarbonyl)-Cyclopento	
	ane
Carboxylic Acid) (A-922500)	ane
Carboxylic Acid) (A-922500)	ane 
Carboxylic Acid) (A-922500) PHOME	ane
Carboxylic Acid) (A-922500)	ane 
Carboxylic Acid) (A-922500)	ane 
Carboxylic Acid) (A-922500)PHOMEPhosphatidylcholine Assay KitPhosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxoyaleryl) Phosphatidylcholine (POV-PC)	ane  nal Antibody)
Carboxylic Acid) (A-922500)PHOMEPhosphatidylcholine Assay KitPhosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxoyaleryl) Phosphatidylcholine (POV-PC)	ane  nal Antibody)
Carboxylic Acid) (A-922500)PHOMEPhosphatidylcholine Assay KitPhosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxoyaleryl) Phosphatidylcholine (POV-PC)	ane  nal Antibody)
Carboxylic Acid) (A-922500)PHOMEPhosphatidylcholine Assay KitPhosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxoyaleryl) Phosphatidylcholine (POV-PC)	ane  nal Antibody)
Carboxylic Acid) (A-922500)	nal Antibody)
Carboxylic Acid) (A-922500) PHOME Phosphatidylcholine Assay Kit Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC) Pinane Thromboxane A <sub>2</sub> PLA <sub>2c</sub> (Adipose Triglyceride Lipase Polyclonal Antibody) POV-PC PPARa Blocking Peptide	nal Antibody)
Carboxylic Acid) (A-922500) PHOME Phosphatidylcholine Assay Kit Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC) Pinane Thromboxane A <sub>2</sub> PLA <sub>2c</sub> (Adipose Triglyceride Lipase Polyclonal Antibody) POV-PC PPARa Blocking Peptide PPARa LBD (human recombinant)	nal Antibody)
Carboxylic Acid) (A-922500)	nal Antibody)
Carboxylic Acid) (A-922500)	ane nal Antibody)23,
Carboxylic Acid) (A-922500). PHOME	ane nal Antibody) 23,
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  sinant))  See PPAI
Carboxylic Acid) (A-922500) PHOME Phosphatidylcholine Assay Kit Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclot 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC) Pinane Thromboxane A2 PLA2. (Adipose Triglyceride Lipase Polyclonal Antibody) POV-PC PPARa Blocking Peptide PPARa Ligand Binding Domain (PPARa LBD (human recomb PPARa Ligand Binding Domain (PPARa LBD (human recomb PPARa Polyclonal Antibody PPARa Transcription Factor Assay Kit PPARB PPARB Blocking Peptide	nal Antibody)
Carboxylic Acid) (A-922500) PHOME Phosphatidylcholine Assay Kit Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC) Pinane Thromboxane A <sub>2</sub> PLA <sub>2c</sub> (Adipose Triglyceride Lipase Polyclonal Antibody) POV-PC PPARa Blocking Peptide PPARa LBD (human recombinant) PPARa Ligand Binding Domain (PPARa LBD (human recomb PPARa Polyclonal Antibody PPARa Transcription Factor Assay Kit PPARB PPARB Blocking Peptide PPARB Blocking Peptide PPARB Blocking Peptide PPARB LBD (human recombinant)	ane  nal Antibody)  23,  pinant))  See PPAI
Carboxylic Acid) (A-922500). PHOME	ane nal Antibody) 23, binant))
Carboxylic Acid) (A-922500). PHOME	nal Antibody) 23, sinant))
Carboxylic Acid) (A-922500). PHOME	nal Antibody) 23, sinant))
Carboxylic Acid) (A-922500). PHOME Phosphatidylcholine Assay Kit Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC) Pinane Thromboxane A2 PLA2c (Adipose Triglyceride Lipase Polyclonal Antibody) POV-PC PPARa Blocking Peptide PPARa LBD (human recombinant) PPARa Ligand Binding Domain (PPARa LBD (human recomb PPARa Polyclonal Antibody	ane  nal Antibody)  23,  sinant))
Carboxylic Acid) (A-922500). PHOME. Phosphatidylcholine Assay Kit. Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclot 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC)	ane nal Antibody)
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody) 23,  sinant))
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  sinant))
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  sinant))  See PPAI
Carboxylic Acid) (A-922500). PHOME Phosphatidylcholine Assay Kit Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC) Pinane Thromboxane A2 PLA2c (Adipose Triglyceride Lipase Polyclonal Antibody) POV-PC PPARa Blocking Peptide PPARa LBD (human recombinant) PPARa Ligand Binding Domain (PPARa LBD (human recomber PPARa Polyclonal Antibody) PPARa Transcription Factor Assay Kit PPARB PPARB Blocking Peptide PPARB Blocking Peptide PPARB Nelsol (human recombinant) PPARB Western Ready Control PPARB Western Ready Control PPARP Blocking Peptide PPARP Blocking Peptide PPARP FL (human recombinant from E. coli) PPARY FL (human recombinant from Sf21 cells) PPARY FUBLEngth	ane  nal Antibody)  23,  sinant))  See PPAI
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  sinant))  See PPAI
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  sinant))  See PPAI
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  pinant))  See PPAI See PPAI  nant))
Carboxylic Acid) (A-922500). PHOME Phosphatidylcholine Assay Kit Phosphatidylcholine-Sterol O-Acyltransferase (LCAT Polyclor 2-(5-oxovaleryl) Phosphatidylcholine (POV-PC) Pinane Thromboxane A2 PLA2, (Adipose Triglyceride Lipase Polyclonal Antibody) POV-PC PPARa Blocking Peptide PPARa Ligand Binding Domain (PPARa LBD (human recombinant) PPARa Polyclonal Antibody PPARa Transcription Factor Assay Kit PPARB PPARB Blocking Peptide PPARB Blocking Peptide PPARB LBD (human recombinant) PPARB Western Ready Control PPARB Western Ready Control PPARB Transcription Factor Assay Kit PPARP FL (human recombinant from E. coli) PPARY FL (human recombinant from \$f21 cells) PPARY FUB Length PPARY LIgand Binding Domain (PPARY LBD (human recombinant) PPARY FUB Length PPARY LBD (human recombinant) PPARY LBD (human recombinant) PPARY LBD (human recombinant) PPARY Transcription Factor Assay Kit	ane  nal Antibody)  23,  pinant))  See PPAI  See PPAI
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  sinant))  See PPAI  mant))  14,
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  pinant))  See PPAI  nant))  See PCSI
Carboxylic Acid) (A-922500). PHOME	ane  anal Antibody)  23,  binant))  See PPAI  anant))  See PPAI  14,  See PCSI
Carboxylic Acid) (A-922500). PHOME	ane  anal Antibody)  23,  binant))  See PPAI  nant))  14,  See PCSI  See IP Recept
Carboxylic Acid) (A-922500). PHOME	ane
Carboxylic Acid) (A-922500). PHOME	ane  anal Antibody)  23,  binant))  See PPAI  See PPAI  nant))  See PCSI  See IP Recept
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  binant))  See PPAI   See PPAI   14,  See PCSI
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)  23,  binant))  See PPAI   See PPAI   14,  See PCSI
Carboxylic Acid) (A-922500). PHOME	ane  nal Antibody)
Carboxylic Acid) (A-922500). PHOME	ane
Carboxylic Acid) (A-922500). PHOME	ane  anal Antibody)  23,  binant))  See PPAI  See PPAI  nant))  See PCSI  See IP Recept
Carboxylic Acid) (A-922500). PHOME	ane  anal Antibody)  23,  binant))  See PPAI  See PPAI  nant))  See PCSI  See IP Recept

6-keto Prostaglandin $F_{1\alpha}$ -d <sub>4</sub> Lipid Maps MS Standard	.38
6-keto Prostaglandin F, ElA Kit	.38
6-keto Prostaglandin F <sub>1α</sub> Quant-PAK	.38
6-keto Prostaglandin F $_{1\alpha}^{1}$ Quant-PAK	.38
Prostaglandin H Synthase 1 (COX-1 Monoclonal Antibody (Clone CX111)).	.16
Prostaglandin H Synthase 2 (COX-2 Polyclonal Antibody (Clone CX229))	
Prostaglandin I Blocking Peptide	
Prostaglandin I Synthase Monoclonal Antibody (Clone isn-1)	.38
Prostaglandin I Synthase (murine) Blocking Peptide	.39
Prostaglandin I Synthase (murine) Polyclonal Antibody	.39
Prostaglandin I Synthase Polyclonal Antibody	.39
Prostaglandin I <sub>2</sub> ReceptorSee IP Recep	tor
Prostaglandin I <sub>2</sub> (sodium salt)	,39
$\begin{array}{lll} & \text{Prostaglandin I}_3^{'} \text{ (sodium salt)} & \\ & \text{15-deoxy-} \Delta^{12,14} \text{-Prostglandin J}_2 & \\ & & \end{array}$	.39
15-deoxy-Δ'-2' -Prostgianain J <sub>2</sub>	.34
PTA <sub>2</sub> (Pinane Thromboxane A <sub>2</sub> )	,33
Renin Fluorogenic Substrate	
Renin Inhibitor Screening Assay Kit	
Resorcinonaphthalein	
trans-Resveratrol	
Rheocyclan (Taprostene (sodium salt))	
RMI 14514 (TOFA)	
Ro 48-8071	
RO1138452 (CAY10441)	
Rosiglitazone	
SAHA	
Salermide	
Scriptaid	
Selexipag (NS-304)	.30
sGC α subunit (Guanylate Cyclase α subunit (soluble) Polyclonal Antibody)	
Polyclonal Antibody)	.18
sGC B subunit (Guanylate Cyclase B subunit (soluble)	
Polyclonal Antibody)	.18
Simvastatin	,4C
Sinthrome® (Acenocoumarol)	
Sintron® (Acenocoumarol)	6
SIRT1 Direct Fluorescence Screening Assay Kit	.27
SIRT1 FRET-Based Screening Assay Kit	
SIRT1 (human recombinant)	.27
SIRT2 Direct Fluorescence Screening Assay Kit	
SIRT2 (human recombinant)	
SIRT3 Direct Fluorescence Screening Assay Kit	
SIRT3 (human recombinant)	
SIRT4 (human recombinant)	
SIRTS (human recombinant)	.27
SIRT6 Direct Fluorescence Screening Assay Kit	
SIRT6 (human recombinant)	
SIRT7 (human recombinant)	
SIRT7 Polyclonal Antibody	.27
SM (Springomyelin Assay Kir)	.42
SMase	
Sodium Butyrate	/2.
Soluble Epoxide Hydrolase Glocking Pepilde	
Soluble Epoxide Hydrolase (FL) Polyclonal Antibody	<del>،4</del> 0 11
Soluble Epoxide Hydrolase (human recombinant)	
Soluble Epoxide Hydrolase Inhibitor Screening Assay Kit	
Soluble Epoxide Hydrolase (murine recombinant from Sf21 cells)	۱۳. ۱۱
Soluble Epoxide Hydrolase Polyclonal Antibody	. <del>-</del> 11
Sphingomyelin Assay Kit	
Sphingomyelinase Assay Kit	
Sphingomyelinase Inhibitor Screening Assay Kit	42
SQ 29,54821,23,	42
SREBF-2	
SREBP-2 Blocking Peptide	
SREBP-2 Cell-Based Translocation Assay Kit	
SREBP-2 Polyclonal Antibody	.42
SREBP-2 Transcription Factor Assay Kit	.43
SREBP-2 Western Ready Control	.42
SRT 1720	
Statin I (Mevastatin)	.26
Stearidonic Acid	
Sterol O-Acyltransferase 1 (ACAT-1)	6
Sterol O-Acyltransferase 2 (ACAT-2)	
Sterol Regulatory Element-Binding Transcription FactorSee SREBI	
T0070907	.34
T0901317	
Taprostene (sodium salt)	,43
3-TDA (3-Thiatetradecanoic Acid)	
Tenovin-1	
Tenovin-6	
5-(Tetradecyloxy)-2-Furoic Acid (TOFA)	
3-Thiatetradecanoic Acid	
2-thio PAF	
THIOTHDOSPOHUIT RECEPTOL	JO.

hromboxane ${\sf A_2}$ Receptor	Receptor
hromboxane B <sub>2</sub>	43
hromboxane B <sub>2</sub> -d <sub>4</sub>	43
hromboxane B <sub>2</sub> 11-dehydrogenase Polyclonal Antiserum	43
hromboxane B <sub>2</sub> EIA Kit	43
hromboxane B2 Express EIA Kit - Monoclonal	43
hromboxane B2 Lipid Maps MS Standard	43
hromboxane B <sub>2</sub> Quant-PAK	
1-dehydro Thromboxane B <sub>2</sub>	44
4411-dehydro Thromboxane B <sub>2</sub> -d <sub>4</sub>	44
1-dehydro Thromboxane B <sub>2</sub> EIA Kit	44
1-dehydro Thromboxane B <sub>2</sub> Lipid Maps MS Standard	44
1-dehydro Thromboxane B <sub>2</sub> Quant-PAK	44
2,3-dinor Thromboxane B <sub>2</sub>	44
2,3-dinor Thromboxane B <sub>2</sub> EIA Kit	44
hromboxane B <sub>3</sub>	44
1-dehydro Thromboxane B <sub>3</sub>	44
hromboxane Synthase Blocking Peptide	44
hromboxane Synthase Polyclonal Antibody	44
IP60 Polyclonal Antibody	29
OFA	
P Receptor (human) Blocking Peptide	
P Receptor (human) Polyclonal Antibody	45
P Receptor (human) Polyclonal FITC Antibody	45
P Receptor (murine) Blocking Peptide	45
P Receptor (murine) Polyclonal Antibody	45
reprostinil	5,45
richostatin A	
riglyceride Assay KitRK 100 (Beraprost (sodium salt))	45
roalitazoneroalitazone	9
ubastatin A (trifluoroacetate salt)	29
1-keto TXB <sub>2</sub> (11-dehydro Thromboxane B <sub>2</sub> ) 1-44069	44
1 -1AD <sub>2</sub> (INFORMOXANE D <sub>3</sub> )	44
J-44069 J-46619	23,45
<sup>17</sup> -U-46619	23,45
J-51605	
J-18666A	
J-61431F (Ciprostene (calcium salt))	1
J-63557A (Furegrelate (sodium salt))	10
Ilmenide (Chenodeoxycholic Acid)	1.4
Imbellatine (Berberine)	14
/asoactive Eicosanoid HPLC Mixture	7
asopressin (Arginine Vasopressin EIA Kit)	40
VARF 42 ((±)-Warfarin)	
±)-Warfarin	40
±)-Warfarin+)-Warfarin	40 11
+)-Warfarin	40 14
Vy 14643	21
K 96480 (Cicaprost)	
oocoumarin ((±)-Warfarin)	۱۵
000001110111 ((±)-111011111	40

# **Item Number Index**

0003	46	19995	-1
0026	34	34004	3
0031		35200	
0044		35210	
0155		35280	
0162	5,45	35370	4
0173	40	38401	1
0174	40	38406	1
0175		38411	
0218		38601	
0240	32	38606	1
0262	27	38611	1
0263		39620	
0264		48001	
0279	30	50210	
0315	27	50211	4
0316		50350	4
0317		50351	
0318		50510	
0341	25	50511	4
0385	27	50650	4
0386		50651	
0495		50861	
0572	27	52400	4
0559	27	52450	4
0582		60402	
0664		60880	
0671	27	60900	3
0699		60902	
3009		60906	
3042		60909	
3079	7	60910	3
3082		60912	
3085		60916	
3086	27	609249,31,	
3121	27	60925	3
3146		60928	
3172		60929	
3174	27	60945	3
3178	27	61700	3
3187		62237	
3220		62935	
3221	14	62945	2
3277	27	70260	
3282	34	70440	1
3284		70510	
3295	27	7051521,	3
	27		3
3295 3452	27 34	7051521, 7054018,	3
3295 3452 3477	27 34 27	70515	2 2
3295 3452 3477 3491	27 34 27 27	70515	3 2 2 3
3295		70515	3 2 3 3
3295 3452 3477 3491		70515	3 2 3 3
3295		70515	3 2 3 3 3
3295		70515	3 2 3 3 3 3
3295		70515	3 2 3 3 3 3
3295		70515	3 2 3 3 3 3 1
3295 3452 3477 3491 3493 3494 3499 3500 3500 3523	27 34 27 27 27 27 27 27 27 27 27 27 27 27 7	70515	3 2 3 3 3 3 1
3295	27 34 27 27 27 27 27 27 27 27 27 27 27 27 7	70515	3 2 2 3 3 3 3 1 4
3295 3452 3477 3491 3493 3494 3499 3500 3504 3523 3523		70515	3223333141
3295 3452 3457 3471 3491 3493 3494 3500 3504 3523 3524 3523 3526 3531		70515	32233331412
3295		70515	322333314122
3295 3452 3457 3471 3491 3493 3494 3500 3504 3523 3524 3523 3526 3531		70515	322333314122
3295		70515	32233333141222
3295 3452 3477 3491 3493 3494 3500 3504 3523 3524 3531 3660 3631		70515	32233333141221
3295 3452 3452 3477 3491 3493 3494 3500 3504 3523 3524 3523 3526 3531 3640 3633 3644		70515	3223333314122211
3295 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3524 3531 3526 3531 3566 3631 3644 3657	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	3 2 2 3 3 3 3 3 1 4 1 2 2 1 1 1 1
3295 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3526 3531 3544 3631 3644 3644 3644 3657		70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 4
3295. 3452. 3452. 3477. 3491 3493. 3494. 3499. 3500. 3504. 3523. 3524. 3531. 3566. 3531. 3644.		70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 4
3295 3452 3452 3477 3491 3493 3494 3500 3504 3523 3523 3526 3531 3560 3631 3644 3657 3686 3789		70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 4 1
3295 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3540 3556 3631 3644 3657 3686 3789	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71800. 71810. 75780. 89430. 89740. 90011. 90110. 90110. 90114. 90165. 90170.	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 4 1 1
3295 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3546 3631 3644 3657 3686 3789 55120	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	3 2 2 3 3 3 3 3 1 4 1 1 1 1 1 1 1 1 1 1 1 1 1
3295 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3523 3526 3531 3666 3631 3644 3657 3686 3789 5120 5210		70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 1 1 3
3295 3452 3452 3457 3491 3493 3494 3499 3500 3504 3523 3526 3531 3546 3531 3566 3631 3644 3657 3686 3789 5120 5210 6440	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	322333331412221111411132
3295 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3523 3526 3531 3666 3631 3644 3657 3686 3789 5120 5210	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	322333331412221111411132
3295 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3540 3566 3631 3644 3657 3686 3789 5120 5210 6440 6440 6440	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 75780. 89430. 89740. 90011. 90114. 901165. 90170. 90210. 90265. 30, 90290. 90310. 18,	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 1 3 2 1
3295 3452 3452 3457 3491 3493 3494 3499 3500 3504 33523 3526 3531 3526 3531 3560 3566 3631 3644 3649 3657 3686 3789 55120 55210 6440 6450 6440	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	32233333141222111141113214
3295 3452 3452 3457 3491 3493 3494 3499 3500 3506 3523 3526 3526 3531 3560 3566 3631 3644 3657 3686 3789 5120 5210 6440 6450 6440 6450 6440 6445	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 1 1 3 2 1 4 1
3295 3452 3452 3457 3491 3493 3494 3494 3499 3500 3504 3523 3526 3531 3566 3631 3644 3657 3686 3789 5120 5210 6440 6445 6445 6465 6460 6465 6465 6481	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 75780. 89430. 89430. 89740. 90011. 90110. 90114. 901165. 90170. 90210. 90265. 30, 90290. 90314. 90320. 90500.	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 1 4 1 1 1 3 2 1 4 1 4
3295 3452 3452 3457 3491 3493 3494 3499 3500 3504 33523 3526 3531 3526 3531 3560 3566 3631 3644 3649 3657 3789 55100 55100 6440 6450 6465 6831 6831	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 4 1 1 1 3 2 1 4 1 4 1
3295 3452 3452 3452 3477 3491 3493 3494 3499 3500 3506 3523 3526 3526 3531 3560 3566 3631 3644 3657 3686 3789 5120 5210 6440 6450 6440 6450 6460 6465 6831 6833 88210	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 1 4 1 1 1 3 2 1 4 1 4 1 3
3295 3452 3452 3457 3491 3493 3494 3499 3500 3504 33523 3526 3531 3526 3531 3560 3566 3631 3644 3649 3657 3789 55100 55100 6440 6450 6465 6831 6831	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	3 2 2 3 3 3 3 3 1 4 1 2 2 2 1 1 1 1 4 1 1 1 3 2 1 4 1 4 1 3
3295 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3566 3531 3566 3631 3644 3657 3886 3789 5120 5210 6440 6440 6445 6465 6460 6465 6831 6833 8210 8211	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70575	32233331412221111411321413
3295 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3526 3531 3540 3566 3631 3644 3657 3686 3789 55120 6440 6445 6456 6450 6450 6450 6450 645	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 75780. 89430. 89740. 90011. 90110. 90110. 90110. 90110. 90110. 90114. 90165. 90170. 90210. 90265. 30, 90290. 90314. 90320. 90500. 100011. 100023. 100027. 100028.	3 2 2 3 3 3 3 3 3 1 4 1 1 2 2 2 1 1 1 1 1 1 1 1 1 3 1 3 1 3 1
3295 3452 3452 3452 3477 3491 3493 3493 3494 3500 3504 3526 3531 3526 3531 3560 3566 3631 3644 3644 3657 3686 3789 55210 6440 6450 6460 6465 6831 6833 8210 8211 8212	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71800. 71810. 75780. 89430. 89730. 89740. 90011. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90114. 90165. 90170. 90265. 30, 90290. 90310. 90320. 90500. 100011 100023. 1000027. 1000028. 1000020.	3223333314122211114111321411331
3295 3452 3452 3452 3457 3491 3493 3494 3499 3500 3504 3523 3526 3531 3566 3631 3666 3631 3644 3657 3886 3789 5120 5210 6440 6445 6465 6460 6465 6831 6833 8210 8211 8212 8213 8215 8215	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71800. 71810. 75780. 89430. 89430. 89740. 90011. 90110. 9	322333314122211114111321413311
3295 3452 3452 3452 3477 3491 3493 3493 3494 3500 3504 3526 3531 3526 3531 3560 3566 3631 3644 3644 3657 3686 3789 55210 6440 6450 6460 6465 6831 6833 8210 8211 8212	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71800. 71810. 75780. 89430. 89730. 89740. 90011. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90114. 90165. 90170. 90265. 30, 90290. 90310. 90320. 90500. 100011 100023. 1000027. 1000028. 1000020.	322333314122211114111321413311
3295 3452 3452 3452 3457 3491 3493 3494 3499 3500 3504 3523 3526 3531 3566 3631 3644 3657 3686 3789 5120 5210 6440 6445 6465 6481 6831 8210 8211 8211 8215 8216 8220	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 75780. 89430. 89430. 89740. 90011. 90110. 90110. 90110. 90110. 90110. 90114. 90165. 90170. 90290. 90314. 90320. 90314. 90320. 90310. 90314. 90320. 90500. 100011. 100023. 100027. 100028. 100003. 10000	32233333141222111141113214113113
3295 3452 3452 3452 3457 3491 3493 3494 3494 3499 3500 3504 3523 3526 3531 3526 3531 3560 3566 3631 3644 3640 6465 6831 6440 6465 6831 8210 8211 8211 8215 8216 8216 8220	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89730. 89740. 90011. 90110. 9	32233331412221114111321414131133
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3560 3566 3631 3644 3657 3686 3631 3646 3657 3686 3631 3646 3631 3646 3631 3789 5120 5210 6460 6465 6831 6833 8210 8211 8212 8213 8216 8220 8220 8230	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71800. 71810. 75780. 89430. 89740. 90011. 90110. 9	3223333314122211114111321413333
3295 3452 3452 3452 3452 3477 3491 3493 3494 3494 3499 3500 3523 3526 3531 3566 3531 3566 3631 3644 3657 3686 3789 5120 6440 6445 6481 6831 8210 8211 8211 8211 8211 8212 8213 8215 8216 8220 8230 8330 8570	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89430. 89740. 90011. 90110. 90110. 90110. 90110. 90110. 90114. 90165. 90170. 90210. 90265. 30, 90290. 90314. 90320. 90314. 90320. 90314. 90320. 90314. 90320. 90000. 100011. 100023. 100027. 100028. 100007. 101700. 33, 101710. 101707. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101707. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700.	322333331412221111411132141131133333
3295 3452 3452 3452 3457 3491 3493 3494 3499 3500 3504 3523 3526 3531 3526 3531 3560 3566 3631 3644 3667 3686 3789 5120 5210 6440 6465 66831 66833 8210 8211 8211 8215 8216 8216 8216 8216 8218 8211 8215 8216 8220 8230 8300 88570 9010	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89730. 89740. 90011. 90110. 9	32233333141222111141113214141311333334
3295 3452 3452 3452 3457 3491 3493 3494 3499 3500 3504 3523 3526 3531 3526 3531 3560 3566 3631 3644 3667 3686 3789 5120 5210 6440 6465 66831 66833 8210 8211 8211 8215 8216 8216 8216 8216 8218 8211 8215 8216 8220 8230 8300 88570 9010	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89430. 89740. 90011. 90110. 90110. 90110. 90110. 90110. 90114. 90165. 90170. 90210. 90265. 30, 90290. 90314. 90320. 90314. 90320. 90314. 90320. 90314. 90320. 90000. 100011. 100023. 100027. 100028. 100007. 101700. 33, 101710. 101707. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101707. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700. 33, 101710. 101700.	32233333141222111141113214141311333334
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3560 35366 3631 3644 3657 3686 3631 3646 3631 3646 3631 3789 5120 5210 6440 6450 6465 6831 8210 8211 8212 8213 8216 8220 8230 8300 8870 9010	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89740. 90011. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 90110. 1010. 90110. 1010. 901	322333314122211141113214141311333342
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3566 3531 3566 3631 3644 3657 3686 3789 5120 5210 6440 6445 6465 6460 6465 6831 8210 8211 8212 8213 8211 8212 8213 8216 8220 8230 8330 8570 9010	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70575	322333331412221111411132141413113333421
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3526 3531 3566 3631 3644 3657 3789 5120 5210 6440 6445 6456 6831 6833 8210 8211 8215 8212 8213 8215 8216 8220 8230 8300 8870 9010	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89430. 89740. 90011. 90110. 9	322333331412221111411132141131133334211
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3526 3531 3560 3566 3631 3644 3657 3686 3631 3644 3657 3686 3631 3646 3631 3646 3631 3886 3789 5120 5210 6440 6455 6450 6465 6831 8210 8211 8212 8213 8216 8210 8211 8212 8213 8216 8220 8230 8300 88570 9010 9020 9021 9020	27 34 27 27 27 27 27 27 27 27 27 27 27 27 34 46 46 41 46 42 27 88 318 38 38 23,45 45 45 5,15 5,10,23 10 10 5,22,23 15,23 5,23,39 5,23,39 5,23,39 3,39 3,34 10,23 23,33 23,33 23,33 23,33 23,33 23,33 21,23,42	70515	322333314122211141113214141311333342111
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3531 3566 3631 3666 3631 3644 3657 3886 3789 5120 5210 6440 6445 6465 6460 6465 6831 8210 8211 8212 8213 8211 8212 8213 8216 8220 8230 8230 8230 8230 8270 9010 9021	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71800. 71810. 75780. 89430. 89430. 89740. 90011. 90110. 9	322333314122211114111321411311333334211111
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3504 3523 3526 3526 3531 3560 3566 3631 3644 3657 3686 3631 3644 3657 3686 3631 3646 3631 3646 3631 3886 3789 5120 5210 6440 6455 6450 6465 6831 8210 8211 8212 8213 8216 8210 8211 8212 8213 8216 8220 8230 8300 88570 9010 9020 9021 9020	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	322333314122211114111321411311333334211111
3295 3452 3452 3452 3452 3477 3491 3493 3494 3494 3499 3500 3504 3523 3526 3531 3566 3631 3644 3657 3686 3789 5120 5210 6440 6445 6485 6481 6831 8210 8211 8212 8213 8215 8216 8220 8230 8330 83670 9010 9020 9020 9020 9025 9030 9050 9380	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89740. 90011. 90110. 90114. 90116. 90110. 90114. 90116. 90110. 90110. 10170. 90210. 90265. 30, 90290. 90314. 90320. 90314. 90320. 90314. 90320. 90001. 100011. 100023. 100027. 100028. 100007. 10170. 10170. 10170. 10170. 10170. 10170. 33, 101710. 101720. 101882. 160070. 160108. 160109. 160100. 160108. 160109. 160110. 160108. 160109. 160110.	3223333141222111411132141413 11333342111111
3295 3452 3452 3452 3452 3477 3491 3493 3494 3499 3500 3506 3523 3526 3523 3526 3531 3560 3566 3631 3644 3657 3686 3631 3644 3657 3686 3631 3646 3631 3640 6450 6460 6450 6460 6450 6460 6450 6460 6450 647 6480 8210 8211 8212 8212 8216 8220 8233 8300 8570 9010 9020 9021 9020 9021 9020 9021 9020 9021 9020 9021 9020 9021 9020 9021 9020 9021 9020 9021 9020 9021 9020 9021	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515	3223333141222111411132141413 113333421111111111111111111111111111111
3295 3452 3452 3452 3452 3477 3491 3493 3494 3494 3499 3500 3504 3523 3526 3531 3566 3631 3644 3657 3686 3789 5120 5210 6440 6445 6485 6481 6831 8210 8211 8212 8213 8215 8216 8220 8230 8330 83670 9010 9020 9020 9020 9025 9030 9050 9380	27 34 27 27 27 27 27 27 27 27 27 27 27 27 27	70515. 21, 70540. 18, 70675. 70730. 70785. 71730. 71740. 71750. 71810. 71810. 75780. 89430. 89740. 90011. 90110. 90114. 90116. 90110. 90114. 90116. 90110. 90110. 10170. 90210. 90265. 30, 90290. 90314. 90320. 90314. 90320. 90314. 90320. 90001. 100011. 100023. 100027. 100028. 100007. 10170. 10170. 10170. 10170. 10170. 10170. 33, 101710. 101720. 101882. 160070. 160108. 160109. 160100. 160108. 160109. 160110. 160108. 160109. 160110.	3223333141222111411132141413 113333421111111111111111111111111111111

5011316	
	10006438
5011616	1000645211,4
5012016	100064801
5012616	1000648210,4
5030425	100065152
5040225	1000652930,4
5060031	10006594
5060231	10006782
3060331	1000679818,3
3060431	10006830
5063038	1000683243,4
5064039	10006855
5070425	10006906
	10006914
6071544	
5072043	10006915
5088029	10006964
5089518	10006998
5089718	10006999
3815011	10007171
0001111	10007185
)170035	10007186
)170733,34,35	10007219
)171033	10007237
1521038	10007239
1903043	10007252
19500	10007253
6007022	10007255
3010616	10007260
5010716	10007261
5040225	10007262
5060031	10007263
	10007274
5060331	
39	10007451
5071544	100075022
5088029	10007599
5088129	10007601
5089518	10007640
5089718	10007663
5090031	100076652
5090631	100076722
0000125	10007685
1512138	10007686
1521138	1000766
1903143	100077671
1905144	100078172
1950144	10007819
3315117	10007820
339518	10007850
3500129	10007852
	1000787511,4
393017	
394018	10007893
	10007893
894018 894516	10007912
39401       8         39451       6         00090       40	10007912 100079231
89401     8       89451     6       00090     40       00150     27	10007912
39401       8         39451       6         00090       40         00150       27         00160       29	10007912
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29	10007912
39401       8         39451       6         00090       40         00150       27         00160       29	10007912
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29	10007912
89401     8       89451     6       90090     40       90150     27       90160     29       90170     29       90230     27       90280     27	10007912. 10007923. 1 10007927. 1 10007941. 3 10008463. 1 10008492. 2 10008596. 4
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27	10007912 10007923
89401     8       897451     6       900090     40       10150     27       10160     29       10170     29       10230     27       10280     27       10290     27       10310     18	10007912
89401     8       89451     6       90090     40       10150     27       10160     29       10170     29       10230     27       10280     27       10290     27       10310     18       10330     42	10007912. 10007923
89401     8       197451     6       100090     40       10150     27       10160     29       10170     29       10230     27       10280     27       10290     27       10310     18       10330     42       10901     30	10007912 10007923
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       50901     30       000183     10	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008596. 40008597. 10008598. 40008599. 40008699. 40008699. 40008699. 40008690. 40008610.
89401     8       197451     6       100090     40       10150     27       10160     29       10170     29       10230     27       10280     27       10290     27       10310     18       10330     42       10901     30	10007912 10007923
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       50901     30       000183     10	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008596. 40008597. 10008598. 40008599. 40008699. 40008699. 40008699. 40008690. 40008610.
89401     8       89451     6       10099     40       10150     27       10160     29       10170     29       10230     27       10280     27       10290     27       10310     18       10333     42       260901     30       100183     10       100462     38       100336     6	10007912. 10007923. 110007927. 10007941
89401     8       197451     6       00090     40       10150     27       20160     29       20170     29       20230     27       20280     27       20290     27       20310     18       20330     42       20090     30       20310     10       200330     42       20091     30       200183     10       200462     38       2000336     6       200337     6	10007912 10007923
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       003330     42       60901     30       000183     10       000462     38       000337     6       000368     15	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008596. 40008597. 10008598. 40008599. 40008609. 40008610. 10008613. 10008811. 400088846. 400088846.
89401     8       89451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       50901     30       000183     10       000462     38       000337     6       000368     15       0004023     43	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008596. 2010008597. 2010008599. 2010008609. 2010008610. 2010008811. 20008811. 20008878. 2010008878. 2010008878. 2010008878. 2010008878. 2010008878. 2010008878.
89401     8       197451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       003330     42       260901     30       000183     10       000462     38       000336     6       000337     6       000368     15       0004023     43       0004110     45	10007912. 10007923. 110007927. 10007941. 10008463. 10008596. 20008597. 20008598. 210008599. 210008609. 210008610. 210008613. 210008811. 210008814. 210008878. 2210008878. 2310008811. 2310008884. 2310008884. 2310008884. 2310008884. 2310008884. 2310008884. 2310008884. 2310008884. 2310008884.
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       60901     30       000183     10       000462     38       000336     6       000368     15       000410     45       0004110     45       0004111     17	10007912 10007923
89401     8       89451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       50901     30       000183     10       000462     38       000336     6       000337     6       000368     15       5004023     43       0004110     45       0004111     17       004380     31	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008597. 10008598. 10008599. 10008609. 10008613. 10008811. 10008846. 10008884. 10008884. 110009884. 110009884. 110009884. 110009884. 110009884. 110009884. 1100090917. 11,5
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       60901     30       000183     10       000462     38       000336     6       000368     15       000410     45       0004110     45       0004111     17	10007912 10007923
89401     8       89451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       50901     30       000183     10       000462     38       000336     6       000337     6       000368     15       5004023     43       0004110     45       0004111     17       004380     31	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008597. 10008598. 10008599. 10008609. 10008613. 10008811. 10008846. 10008884. 10008884. 110009884. 110009884. 110009884. 110009884. 110009884. 110009884. 1100090917. 11,5
89401     8       897451     6       900090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       003330     42       60901     30       000183     10       000462     38       000336     6       000337     6       000368     15       000410     45       0004110     45       0004452     31       0004110     45       0004452     45       0004452     45       0004466     16	10007912 10007923 10007927 10007921 10007941 10008463 10008492 10008596 20008597 20008598 20008699 20008610 10008610 1000881 10008884 10008888 10008888 10009017 11,10009085 10009088 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009085 10009088 10009088
89401     8       89451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       003330     42       50901     30       000183     10       000462     38       000337     6       000338     15       0004023     43       0004110     45       0004111     17       0004380     31       0004946     16       0004971     16	10007912. 10007923. 110007927. 10007941. 10008463. 10008596. 10008597. 10008598. 10008609. 110008610. 110008613. 10008811. 10008884. 10008884. 110009888. 110009888. 110009888. 110009915. 1100090918. 1100090918. 1100090918. 1100090918. 11000901914. 110009110009134. 110009134. 110009127. 110009134. 110009145. 110009141.
89401     8       89451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00333     42       40901     30       000183     10       000462     38       000336     6       000337     6       000368     15       0004023     43       0004110     45       0004111     17       0004380     31       0004971     16       0005056     17	10007912. 10007923. 110007923. 120007924. 10007941. 10008463. 10008596. 10008597. 10008599. 10008609. 10008610. 110008811. 10008818. 10008878. 10008884. 110008884. 110009883. 120009883. 120009917. 111.31000988. 120009918. 120000918. 12000000000000000000000000000000000000
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       003330     42       60901     30       000183     10       000462     38       000336     6       000337     6       000368     15       000410     45       0004111     17       0004452     45       0004946     16       0004971     16       0005056     17       0005090     6	10007912. 10007923. 110007927. 10007941. 10008463. 10008596. 201008599. 201008609. 201008610. 201008811. 201008811. 201008884. 201008884. 201009098. 201009915. 20100915. 20100915. 2010009231. 201000927. 20100927. 20100915. 20100915. 201000927. 20100915. 201000927. 201000915. 201000927. 201000927. 201000915. 201000927. 201000927. 201000927. 201000915. 201000915. 201000915. 201000927. 2010000927. 20100000000000000000000000000000000000
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       003330     42       50901     30       000183     10       000462     38       000337     6       000368     15       000410     45       000411     17       0004380     31       0004966     16       0004971     16       0005090     6       0005091     6	10007912. 10007923. 110007927. 10007941. 10008463. 10008596. 10008597. 10008598. 10008599. 10008609. 110008610. 110008811. 10008884. 10009878. 10008884. 10009017. 11,510009085. 10009088. 10009018. 10009019. 11,0009019.
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       003330     42       60901     30       000183     10       000462     38       000336     6       000337     6       000368     15       000410     45       0004111     17       0004452     45       0004946     16       0004971     16       0005056     17       0005090     6	10007912. 10007923. 110007927. 10007941. 10008463. 10008596. 201008599. 201008609. 201008610. 201008811. 201008811. 201008884. 201008884. 201009098. 201009915. 20100915. 20100915. 2010009231. 201000927. 20100927. 20100915. 20100915. 201000927. 20100915. 201000927. 201000915. 201000927. 201000927. 201000915. 201000927. 201000927. 201000927. 201000915. 201000915. 201000915. 201000927. 2010000927. 20100000000000000000000000000000000000
39401       8         39451       6         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         50901       30         000183       10         000462       38         000337       6         000337       6         000338       15         0004023       43         0004110       45         0004111       17         0004380       31         0004974       16         0005056       17         0005090       6         0005091       6         0005186       10,23	10007912. 10007923. 110007923. 110007927. 10007941. 10008463. 10008492. 10008596. 201008599. 201008609. 201008610. 20100861. 20100861. 201008681. 201008681. 201008681. 201008681. 201008681. 201008884. 201008884. 201009917. 201008884. 201009918. 201009929. 201009929. 2010009299. 2010009299.
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00333     42       60901     30       000183     10       000336     6       000337     6       000368     15       000410     45       000411     17       004380     31       0004971     16       0005056     17       0005090     6       0005186     10,23       0005186     10,23       0005257     39	10007912. 10007923. 110007927. 10007941. 10008463. 10008596. 20008597. 20008599. 20008609. 20008610. 20008613. 20008811. 20008884. 20008884. 20009088. 20009088. 20009088. 20009098.
89401     8       897451     6       00090     40       00150     27       00160     29       00170     29       00230     27       00280     27       00290     27       00310     18       00330     42       80901     30       000183     10       000462     38       000337     6       000368     15       000410     45       000411     17       000482     45       0004946     16       0005971     16       0005090     6       0005091     6       0005257     39       0005263     45	10007912. 10007923. 110007927. 10007941. 10008463. 10008596. 201008598. 201008599. 201008610. 20100861. 20100861. 201008684. 201008884. 201008884. 201008884. 201009888. 201009888. 201009988. 201009988. 2010099988. 2010099231. 2010099286. 2010099286. 201000927. 201009286. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928. 201009928.
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         50901       30         000183       10         000462       38         000337       6         000338       15         0004023       43         0004110       45         0004111       17         0004380       31         0004974       16         0005971       16         0005091       6         000556       17         0005091       6         0005263       45         0005263       45         0005292       8	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008597. 10008598. 10008599. 10008610. 11000881. 10008818. 10008884. 10008884. 10009017. 11,510009088. 1000909088. 100090909090909090909090909090909090909
39401       8         39451       6         39451       6         000990       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00333       42         50901       30         000183       10         000462       38         000336       6         000337       6         000368       15         0004023       43         0004110       45         0004111       17         0004380       31         0004971       16         0005090       6         0005091       6         0005091       6         0005292       8         0005368       17,34	10007912. 10007923. 10007923. 10007927. 10007941. 33 10008463. 10008596. 20 10008597. 21 10008599. 21 10008609. 22 10008610. 21 10008611. 21 10008811. 21 20008884. 21 20009017. 21 200909088. 22 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016.
39401       8         39451       6         39451       6         000990       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00333       42         50901       30         000183       10         000462       38         000336       6         000337       6         000368       15         0004023       43         0004110       45         0004111       17         0004380       31         0004971       16         0005090       6         0005091       6         0005091       6         0005292       8         0005368       17,34	10007912. 10007923. 110007927. 10007941. 10008463. 10008492. 10008597. 10008598. 10008599. 10008610. 11000881. 10008818. 10008884. 10008884. 10009017. 11,510009088. 1000909088. 100090909090909090909090909090909090909
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         60901       30         000183       10         000462       38         000336       6         000368       15         000410       45         000411       17         0004380       31         0004946       16         0005056       17         0005090       6         0005186       10,23         0005257       39         0005263       45         0005518       17,34         0005518       17,34	10007912. 10007923. 10007923. 10007927. 10007941. 33 10008463. 10008596. 20 10008597. 21 10008599. 21 10008609. 22 10008610. 21 10008611. 21 10008811. 21 20008884. 21 20009017. 21 200909088. 22 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009015. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016. 20 2009016.
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         003330       42         50901       30         000183       10         000462       38         000337       6         000338       15         0004010       45         000411       17         0004380       31         0004452       45         0004946       16         000597       6         0005091       6         0005091       6         0005263       45         0005292       8         0005518       17,34         0005519       22	10007912. 10007923. 10007927. 10007941
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         50901       30         000183       10         000462       38         000336       6         000337       6         0004023       43         0004110       45         0004111       17         0004380       31         0004974       16         0005056       17         0005090       6         000527       39         0005263       45         0005292       8         0005368       17,34         0005519       22         0005745       15,34	10007912. 10007923. 10007927. 10007941. 33 10008463. 10008492. 10008596. 201008599. 201008609. 201008610. 201008613. 201008811. 201008811. 201008878. 201008878. 201008878. 201008884. 201009017. 201008846. 201009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 2010009018. 201000018. 2010000018. 20100000000000000000000000000000000000
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         60901       30         000183       10         000462       38         000336       6         000338       15         0004023       43         0004110       45         0004111       17         0004380       31         0004946       16         0005056       17         0005090       6         0005091       6         0005263       45         0005292       8         0005518       22         0005519       22         0005745       15,34         0005913       10,23	10007912. 10007923. 10007923. 10007927. 10007941. 33 10008463. 10008596. 40008597. 40008599. 40008609. 40008610. 40008613. 40008811. 40008811. 40008884. 40009017. 40009088. 40009018.
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00330       42         50901       30         000183       10         000462       38         000337       6         000338       15         00040403       43         0004110       45         0004111       17         0004380       31         0004946       16         0004971       16         0005090       6         0005091       6         0005292       8         0005293       8         0005297       8         000518       17,34         0005519       22         0005913       10,23         0005913       10,23         0005913       10,23         0005913       10,23         0005913       10,23         0005914       19	10007912. 10007923. 10007927. 10007941. 201007941. 201008596. 201008597. 201008599. 201008610. 20100861. 20100861. 20100861. 20100861. 20100861. 20100861. 20100861. 20100861. 20100861. 20100861. 201008846. 201008846. 201008884. 201009888. 201009915. 2010009018. 201009018. 201009019. 2010009019. 2010009019. 2010009019. 2010009019. 2010009019. 2010000019. 201000019. 2010000
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         50901       30         000183       10         000462       38         000337       6         000337       6         000348       15         0004023       43         0004110       45         0004111       17         0004380       31         0004946       16         0004971       16         000506       17         0005090       6         0005263       45         0005263       45         0005518       10,23         0005519       22         0005745       15,34         0006217       39	10007912. 10007923. 10007927. 10007941. 33 10008463. 10008492. 10008596. 40008597. 40008599. 40008610. 40008610. 40008611. 40008611. 40008681. 40008878. 40008884. 40009017. 40008884. 40009017. 40009018. 400
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         003330       42         60901       30         000183       10         000462       38         000336       6         000337       6         000368       15         000410       45         0004111       17         0004380       31         0004971       16         0005056       17         0005090       6         0005186       10,23         0005257       39         0005263       45         0005718       22         0005518       22         0005518       22         0005745       15.34         0006217       34         0006247       34	10007912. 10007923. 10007923. 10007927. 10007941. 33 10008463. 10008492. 10008596. 20008597. 20008599. 20008609. 2010008610. 2010008613. 201008811. 201008811. 201008884. 201009884. 201009917. 2010098884. 201009918. 20100
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         003330       42         50901       30         000183       10         000462       38         000337       6         000338       15         0004010       45         000411       17         0004380       31         0004452       45         0004946       16         000597       6         0005091       6         0005091       6         0005263       45         0005292       8         0005518       17,34         0005519       22	10007912. 10007923. 10007927. 10007941. 33 10008463. 10008492. 10008596. 40008597. 40008599. 40008610. 40008610. 40008611. 40008611. 40008681. 40008878. 40008884. 40009017. 40008884. 40009017. 40009018. 400
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         003330       42         60901       30         000183       10         000462       38         000336       6         000337       6         000368       15         000410       45         0004111       17         0004380       31         0004971       16         0005056       17         0005090       6         0005186       10,23         0005257       39         0005263       45         0005718       22         0005518       22         0005518       22         0005745       15.34         0006217       34         0006247       34	10007912. 10007923. 10007923. 10007927. 10007941. 33 10008463. 10008492. 10008596. 20008597. 20008599. 20008609. 2010008610. 2010008613. 201008811. 201008811. 201008884. 201009884. 201009917. 2010098884. 201009918. 20100
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         00901       30         000183       10         000462       38         000337       6         000338       15         0004023       43         0004110       45         0004111       17         0004380       31         0004946       16         0004971       16         0005056       17         0005090       6         0005263       45         000527       39         0005263       45         0005186       10,23         0005519       22         0005745       15,34         000627       39         000627       39         000627       39         000627       39         000627	10007912. 10007923. 10007923. 110007927. 10007941. 10008463. 10008492. 10008596. 201008597. 201008599. 201008609. 201008610. 201008610. 201008611. 201008611. 201008611. 201008611. 201008611. 201008611. 201008611. 201008811. 201008811. 201008811. 201008884. 2010008884. 2010009017. 201008884. 201009018. 201009018. 2010009018. 2010009018. 201009018. 201009018. 201009018. 201009018. 201
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         60901       30         000183       10         000462       38         000336       6         000337       6         000368       15         000410       45         004111       17         004380       31         004946       16         0005056       17         0005090       6         0005186       10,23         000527       39         0005263       45         000591       6         000518       17,34         0005918       22         0005745       15,34         0005745       15,34         0005913       10,23         000527       39         0006217       39         000	10007912. 10007923. 10007923. 110007927. 10007941. 10008463. 10008492. 10008596. 201008599. 201008609. 10008610. 201008613. 201008811. 201008811. 201008811. 201008884. 201009884. 201009917. 201009884. 201009918. 2010009918. 201000918. 201
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         50901       30         000183       10         000462       38         000336       6         000368       15         000410       45         000411       17         0004380       31         0004946       16         0005091       6         0005091       6         0005091       6         0005092       8         0005257       39         0005263       45         0005745       15,34         0005913       10,23         000518       17,34         0005913       10,23         0005745       15,34         0005745       15,34         0005745       15,34         0006277       39	10007912. 10007923. 10007927. 10007941. 10008463. 10008596. 201008598. 201008610. 20100861. 20100861. 2010086884. 201008884. 201008884. 201009888. 201009988. 201009888.
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00333       42         50901       30         000183       10         000462       38         000337       6         000338       15         0004010       45         000411       45         000411       17         0004380       31         0004946       16         0004971       16         0005090       6         0005091       6         0005292       8         0005297       8         0005297       8         000518       17,34         000591       6         0005929       8         000519       22         0005745       15,34         000591       22         000591       22         000591	10007912. 10007923. 10007927. 10007941. 2007941. 2008596. 2008596. 2010008597. 2010008599. 2010008610. 201000861. 201008811. 201008884. 2010008884. 2010008884. 2010009017. 201009018. 201009019. 2010009019. 2010009019. 2010009019. 2010009019. 2010009019. 2010009019. 2010009019. 2010009019. 2010009
39401       8         39451       6         00090       40         00150       27         00160       29         00170       29         00230       27         00280       27         00290       27         00310       18         00330       42         60901       30         000183       10         000462       38         000336       6         000337       6         000368       15         000410       45         000411       17         004380       31         004946       16         000556       17         0005090       6         0005186       10,23         0005292       8         0005293       8         0005294       8         0005295       8         0005296       7         0006297       13         000591       6         0005292       8         0005293       8         0005294       10         22       10	10007912. 10007923. 10007927. 10007941. 10008463. 10008596. 201008598. 201008610. 20100861. 20100861. 2010086884. 201008884. 201008884. 201009888. 201009988. 201009888.

1000979727
1000979827
1000987011
1000987132
1000988019,34
1000989311
10009926
10007728
10009929
1000998440
1000998735
1000999540
1000999640
1001008131
1001009616
1001014641
1001014741
100101538
1001022831
1001022931
1001030345
100103327
1001033714,18
10010338
1001033914,28
10010337
10010340
1001034414,40
1001039621,23
1001040127
1001041130
1001041228
1001048616
100105179
1001056727
100105696
1001097128
1001099127
1001102027
1001105423
1001112524
100111319
10011131
1001119127
1001119427
1001121119,34
1001123614
1001128614
1001129618
1001134823,43
1001156221,23,24
1001156327
1001156427
1001156627
1001166940
1001167041
1001167141
1001244714,23
1001244714,23 1001253634
10012447
1001244714,23 1001253634