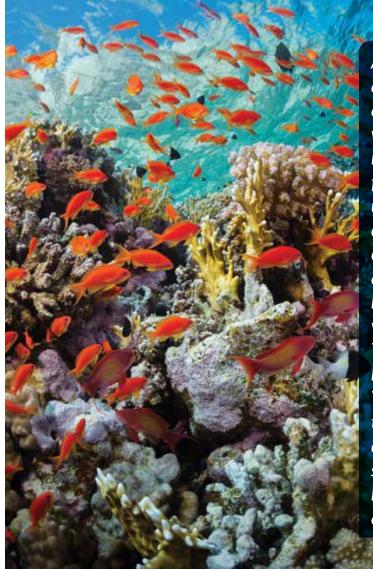


# MEASURE WITH CONFIDENCE

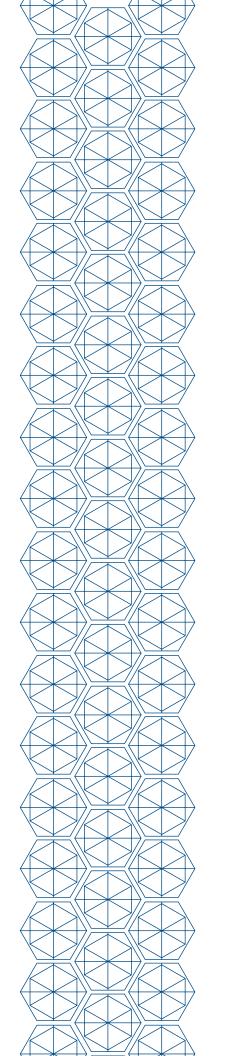


As a direct assay kit manufacturer, we have a unique understanding of assay development and performance. With our assay kits, you will receive all necessary reagents, plates, instructions, and analysis tools to help you find answers quickly and efficiently. We understand that your research is being done with finite time and money, so we create and refine our kits to be affordable and to provide fast, accurate results. You have access to more than 350 assay kits in various formats, including ELISA, cell-based, TR-FRET, enzyme activity, inhibitor screening, high-throughput screening, and transfection. These kits are backed by a team of more than 100 in-house scientists, the same who develop our products and provide technical support, as well as state-of-the-art laboratories, a rigorous validation process, and a wealth of technical data.



# TABLE OF CONTENTS

- 2 Immunoassays to Mass Spectrometry
- **3** Assay Kit Formats
  - Enzyme-Linked Immunosorbent Assays
  - Cell-Based Assays
  - Activity & Detection Assays
  - Related Tools
- **8** Bioanalytical Assay Services
- **9** Kits by Research Area
  - Cancer
  - Cardiovascular Disease
  - Cell Biology
  - Eicosanoids & Lipids
  - Epigenetics
  - Immunology & Inflammation
  - Metabolism
  - Neuroscience
  - Nitric Oxide & Oxidative Injury
  - Toxicology
- **15** Kit Information (listed alphabetically)
- 77 Sample Preparation & Screening Libraries
- **79** Index

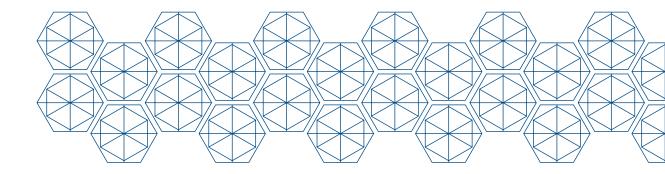


# IMMUNOASSAYS TO MASS SPECTROMETRY: TRANSITIONING TECHNOLOGY

In 1987, there was a need for an easy-to-use, low-waste, and cost-effective technique for the detection and quantification of small bioactive molecules. Cayman Chemical responded to the research community by introducing one of the first commercially-available enzyme-linked immunosorbent assays (ELISAs) for prostaglandin  $\rm E_2$  as an alternative to the cumbersome radioimmunoassay. Almost 30 years later, Cayman continues to offer several ELISAs to researchers, constantly improving on the reliability and sensitivity of these assays.

It is known that antibody-based techniques are limited by the specificity of the antibodies involved, which remains a challenge for researchers quantifying small molecules and biomolecules such as eicosanoids, proteins, and peptides. Moreover, quantitation is limited to the antibody-specific antigen(s), usually a single analyte. Researchers and clinicians alike are now turning to mass spectrometry (MS) as an alternative analytical technique because of its advantages over ELISA-based methods. MS offers an inherent specificity advantage as well as multiplexing capabilities, albeit with significant start-up costs. The value of rapidly quantifying multiple targets simultaneously with accuracy and certainty makes MS an increasingly popular method of choice.

Cayman is committed to leading efforts to develop MS-based approaches, as we have done with ELISAs, and offers a variety of highly pure standards for analysis by MS. Cayman's team of scientists is dedicated to both ELISA and MS analyses to provide researchers worldwide with the tools and expertise needed to make research possible.



#### **ELISAs**

Enzyme-linked immunosorbent assays (ELISAs), also known as enzyme immunoassays (EIAs), combine antibody binding with enzymatic detection to quantify molecules of interest. Cayman Chemical manufactures more than 130 easy-to-use ELISAs in both sandwich/immunometric and competitive formats. Each kit undergoes extensive testing and is backed by a team of in-house scientists committed to offering high-quality products at affordable prices. Many of Cayman's ELISAs employ acetylcholinesterase (AChE), a highly stable enzyme that offers the following advantages over other enzyme reporters:

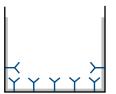
- Does not self-inactivate during turnover
- Enables redevelopment if assay is accidentally splashed or spilled
- Wide pH range (pH 5-10)
- Not affected by buffer salts or preservatives
- A stop reagent is not required
- Plate may be read at any time

#### Cayman's ELISAs also offer:

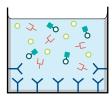
- Precise, sensitive readouts
- Low inter- and intra-assay variability
- Data analysis tools to aid with calculations
- World-class technical support backed by Cayman scientists

Competitive ELISAs are most commonly used to measure small molecules (i.e., lipids, hormones, and small peptides). They are based on the competition between the analyte of interest and an enzyme-conjugated version of the same analyte (tracer) for a limited number of specific antibody binding sites. The amount of tracer that binds to the antibody is inversely proportional to the concentration of the analyte in question.

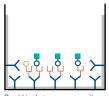
Sandwich/Immunometric ELISAs are suitable for proteins and peptides greater than 20 amino acids in length. This assay format involves two antibodies that "sandwich" the analyte in question between them. A specific capture antibody coats the assay plate, and a second detection antibody is added to the well. The amount of analyte present is directly proportional to the amount of detection antibody, which is typically quantified by its enzyme conjugate.

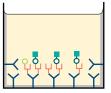


capture antibody and blocked with a proprietary formulation of proteins.



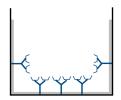
antibody, and either standard or sample



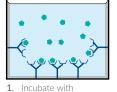




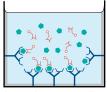
Competitive ELISA



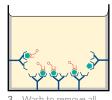
Wells are pre-coated with capture antibody and blocked with a proprietary formulation of proteins.



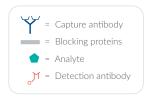
standard or sample



unbound reagents. Incubate with the detection antibody



unbound reagents. stop solution.



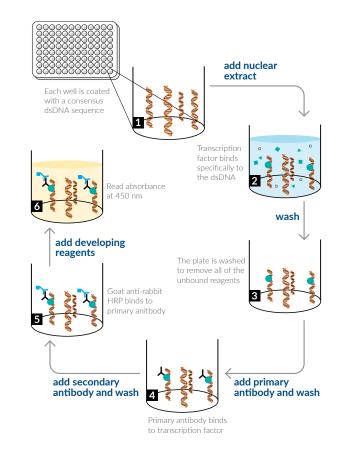
Sandwich/Immunometric ELISA

#### ightarrow Transcription Factor Assays

Cayman Chemical's transcription factor assays are a colorimetric, non-radioactive, and sensitive method for detecting specific transcription factor DNA binding activity in nuclear extracts. This assay format replaces the cumbersome and unreliable electrophoretic mobility shift assays historically used to detect and quantify transcription factors.

This assay format immobilizes a dsDNA sequence containing the target-specific response element onto the wells of a 96-well plate. Transcription factors contained in a nuclear extract bind specifically to the DNA at the response element and are detected by a specific primary antibody. A secondary antibody conjugated to HRP is added to provide a colorimetric readout at 450 nm.

A Nuclear Extraction Kit (Item No. 10009277) is available as a companion to Cayman's transcription factor kits to aid in the isolation of nuclear and cytoplasmic fractions from cultured cells or tissue homogenates. (see page 77)



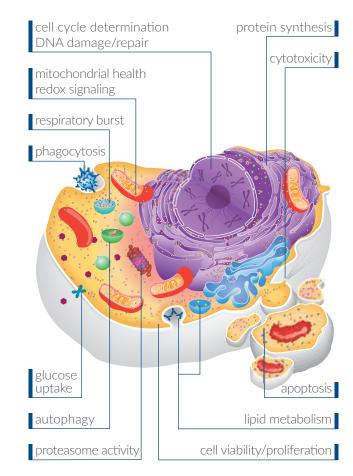
Transcription Factor Assay Protocol

# **CELL-BASED ASSAYS**

Cayman Chemical offers more than 60 cell-based assays for use in studying a variety of cellular functions. Performed in a biologically relevant context, cell-based assays offer a flexible and adaptable method to detect perturbations in function following treatment with a compound of interest. Many of the kits follow simple protocols and can be used with multiple readout formats, making these assays highly adaptable across labs and amenable to high-throughput screening applications.

Kits are available to measure:

- Basic cellular metabolism: uptake of glucose to protein synthesis, oxygen consumption, and glycolysis
- Cell viability/death: cell proliferation, viability, and apoptosis
- Neutrophil and monocyte functions: NET formation, enzyme activity, and phagocytosis
- Lymphocyte function: proliferation and cytotoxicity
- Atherosclerosis and obesity: cellular lipid uptake, trafficking and synthesis, and adipocyte formation



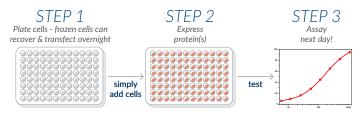
Assays Available for Studying a Variety of Cellular Functions

#### **Reverse Transfection Assays**

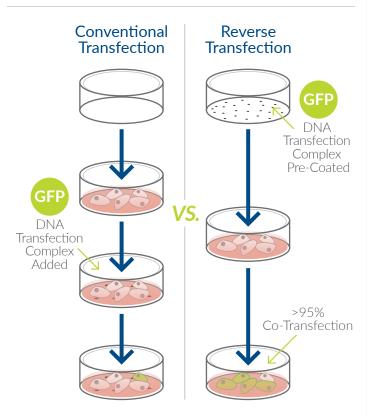
Reliable, reproducible transfection assays in a format that's both convenient and scalable.

- No need to develop, maintain, or expand stable cell lines
- Avoid cell lineage variability
- Suitable for difficult targets for which stable cell lines are not sustainable

Reverse Transfection Reporter Assay technology uses plates pre-coated with optimized transfection complexes, offering up to 3-fold increased transfection and >95% co-transfection efficiency that's consistent well to well. This unique system is ideal for rapid assay development and high-throughput screening.



Reverse Transfection Assay Setup



Comparison of Transfection Methods

#### **ACTIVITY & DETECTION ASSAYS**

Cayman Chemical's activity and detection assays are designed to quantify biomarkers for biological processes, assess mitochondrial function/toxicity, or screen for compounds that affect the activity of various enzymes.

# **Detection Assays**

Cayman Chemical offers a robust line of detection assays for use in metabolism, oxidative stress, immunology, and epigenetics research. These assays have been developed for quick and accurate quantification of biomarkers including small molecules, proteins, and signaling ions such as calcium and phosphate. These assays follow easy, two- to three-step protocols consisting of short incubations with colorimetric or fluorometric substrates followed by a kinetic or end-point analysis. Kits have been developed for use on standard 96-well plate readers and include detailed instructions, all necessary reagents, and sample preparation protocols to run your analysis.

- Simple and fast protocols
- Accurate and reproducible
- Low intra- and inter-assay variability
- Validated by in-house Cayman scientists for relevant matrices
- Backed by world-class technical support

# Activity and Compound Screening Assays

Cayman Chemical offers more than 50 unique in vitro assays to screen enzyme activity. Optimized to save you time in assay development, these assays are easy to use, sensitive, and reliable options to help researchers evaluate the effect of compounds on target activity. Kits come equipped with all the necessary components including detailed protocols, active enzymes, substrates, and specific inhibitors to allow you to run assays with speed and confidence.

#### MitoCheck® Bioenergetics Assays

Create a full mitochondrial function and toxicity profile with simple and cost-effective MitoCheck® assay kits.

- No need for expensive equipment or lengthy incubations
- Adapted for use with microplate readers
- Amenable to high-throughput screening

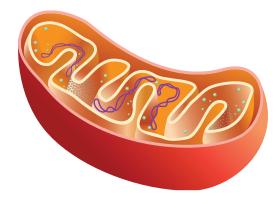
MitoCheck® kits offer researchers the simplest and most biochemically relevant system for studying mitochondrial function. Designed as a set of tools to evaluate mitochondrial toxicity, the kits can function together or as stand-alone assays. Assays are available to screen potential inhibitors of the electron transport chain, assess mitochondrial function in cells by detecting the rate of oxygen consumption, determine citrate synthase activity, and measure mitochondrial membrane potential.

To learn more about methods for mechanism-based mitochondrial toxicity screening, read our application note: Screening of Antipsychotic Compounds for Mitochondrial Toxicity. (www.caymanchem.com/mitonote)

# **Primary Screening Assays** (OCR, membrane potential)



(Complexes I-V)



MitoCheck® Assays

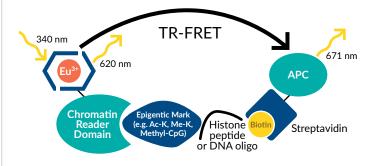


Cayman Chemical's TR-FRET assays are designed to characterize the interaction of chromatin reader domains with potential inhibitors or probes.

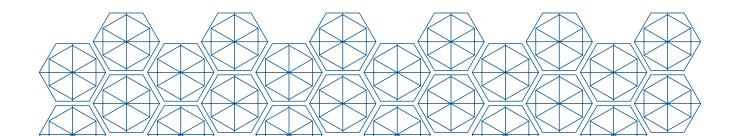
- Rapid and amenable to high-throughput screening
- Homogeneous mix and read TR-FRET in 384-well format
- Substrate-independent, non-kinetic assay
- Miniaturized to final assay volume of 20 μl

Fluorescence Resonance Energy Transfer (FRET), a useful method for detecting protein/protein interactions, is based on the transfer of energy between distinct donor and acceptor fluorophores. When in close proximity (~100 Å), excitation of the donor fluorophore triggers an energy transfer to the acceptor fluorophore, which then emits fluorescence at a distinct wavelength. TR-FRET employs the principles of FRET, but uses a donor fluorophore with a significantly longer fluorescent half-life. This allows for a time delay to be introduced between excitation of the donor fluorophore and the fluorescence measurement. This method minimizes background fluorescence interference, which is particularly useful for high-throughput screening.

In Cayman's TR-FRET assays, the donor fluorophore is a specific recombinant reader domain labeled with a europium chelate. A biotinylated peptide containing a target acetylated lysine serves as a ligand for the reader domain. APC-labeled streptavidin binds with high affinity to the peptide substrate via the biotin moiety and serves as the acceptor fluorophore. Inhibition of the bromodomain/peptide interaction results in loss of the TR-FRET signal.



TR-FRET Assay

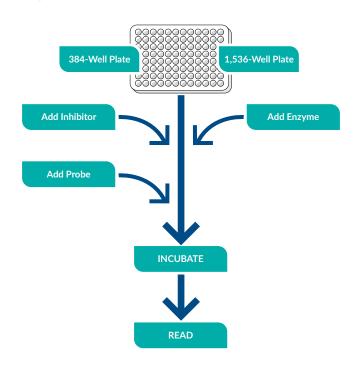


#### SAM-Screener™ Assays

Screen for SAM-competitive methyltransferase inhibitors with SAM-Screener $^{\text{TM}}$  one-step, fluorescent polarization assays.

- Non-kinetic, substrate-independent, and scalable to 1,536-well format
- Plate-based fluorescence polarization measurement
- Assay kits currently available for SET7/9, GLP, and MLL1 enzymes
- Amenable to high-throughput screening

S-Adenosyl-L-methionine (SAM) is a co-substrate of lysine methyltransferases utilized during the methylation of lysine residues on various proteins, including histones. Cayman Chemical's SAM-Screener<sup>™</sup> assays provide a proprietary small molecule fluorescent probe that binds in the SAM binding pocket. Binding of the probe induces a pronounced increase in fluorescence polarization (FP), whereas displacement of the probe by SAM-binding site inhibitors results in concentration-dependent loss of the FP signal. The SAM-Screener<sup>™</sup> assay platform offers a method for studying substrate recognition by methyltransferases and identifying SAM-competitive inhibitors. This quick and efficient assay eliminates the need for complicated multi-step or radioactive enzyme assays, thereby allowing rapid screening of large chemical libraries for new and specific methyltransferase inhibitors.



SAM-Screener™ Assay Setup

# RELATED TOOLS Sample Preparation

Sample preparation is the first and most important step in any assay procedure. Cayman Chemical makes this process quick and convenient by packaging necessary reagents and/or materials plus simplified instructions for the isolation, purification, or extraction of various experimental samples. These products are designed to help circumvent common errors that may occur during the preparation of experimental samples and serve as companions to various assay kits included in this catalog.

For a complete list, see page 77.

#### ightarrow Screening Libraries

Cayman Chemical offers a variety of chemical libraries produced from our collection of high-quality and high-purity compounds. Designed for use in assay screening and drug discovery, focused plates are available for bio-active lipids, fatty acids, prostaglandins, kinase inhibitors, epigenetic compounds, and stem cell small molecule screening applications. The screening libraries are available in a 96-well Matrix tube rack format as stock solutions in DMSO. Any potential screening hit may be purchased individually for activity confirmation and further analysis.

Custom libraries produced from Cayman's compound collection of more than 6,000 molecules are also available to meet your exact screening requirements.

For a complete list, see page 78.



#### **Technology Partners**

Cayman Chemical is proud to partner with the following companies in delivering high-quality research products:











# CONTRACT RESEARCH SERVICES FROM CAYMAN



We offer you the deep and diverse knowledge of more than 100 in-house scientists and the expertise that comes from more than three decades of designing and developing the tools to advance biomedical research.

Cayman Chemical is known as a partner to scientists worldwide, responding to your needs with services in:

- Analytical chemistry
- Assay development
- Bioanalytical assay
- Bioenergetics screening
- Custom antibody services
- Custom organic synthesis
- Epigenetic screening and profiling
- Medicinal chemistry
- Protein production and purification
- Structural biology

# BIOANALYTICAL ASSAY SERVICES



Don't have a plate reader, time, or the resources available to run your assays?

Cayman Chemical's bioanalytical assay services support your needs for drug discovery and development, biomarkers, and basic research. We offer complete sample analysis for analytes measured by our off-the-shelf assays. With our line of well-characterized immunoassays, cell-based assays, and biochemical kits, combined with our knowledgeable and highly experienced scientific staff, we provide accurate, high-quality data. Our bioanalytical assay services include new assay development, customized method development, and sample analysis in multiple biological matrices. As a direct assay kit manufacturer, our staff has a unique understanding of assay development and performance.

To learn more about these services, or to discuss your contract research service needs, please contact us.

Email us at contractresearch@caymanchem.com Visit us online at www.caymanchem.com/services Call us toll free at (888) 526-5351

# KITS BY RESEARCH AREA



7-AAD/CFSE Cell-Mediated Cytotoxicity Assay Kit	1
Acid Phosphatase Colorimetric Assay Kit	1
Aldehyde Dehydrogenase Activity Assay Kit	1
Annexin V FITC Assay Kit	1
Antioxidant Assay Kit	1
Ascorbate Assay Kit	18
Autotaxin Inhibitor Screening Assay Kit	1
Caspase-3 Fluorescence Assay Kit	2
CFSE Cell Division Assay Kit	2
COX Activity Assay Kit	2
COX Colorimetric Inhibitor Screening Assay Kit	2
COX (ovine) Colorimetric Inhibitor Screening Assay Kit	2
COX Fluorescent Activity Assay Kit	
COX Fluorescent Inhibitor Screening Assay Kit	
COX (ovine) Inhibitor Screening Assay Kit	2
COX (ovine/human) Inhibitor Screening Assay Kit	
COX-1 (human) Inhibitor Screening Assay Kit	
COX-1 (ovine) Inhibitor Screening Assay Kit	
COX-2 (human) Inhibitor Screening Assay Kit	
COX-2 (ovine) Inhibitor Screening Assay Kit	
DNA Methylation ELISA Kit	
DNA/RNA Oxidative Damage ELISA Kit	2
DPP (IV) Inhibitor Screening Assay Kit	
EP <sub>4</sub> Receptor (rat) Activation Assay Kit (cAMP)	
EP <sub>4</sub> Receptor (rat) Reporter Assay Kit	
GLP SAM-Screener™ Assay Kit	
Glucose Uptake Cell-Based Assay Kit	
Glutathione Assay Kit	3
Glutathione Peroxidase Assay Kit	3
Glyco-SPOT DNA Repair Assay Kit	
HAT Inhibitor Screening Assay Kit	
HDAC Cell-Based Activity Assay Kit	3
HDAC Fluorometric Activity Assay Kit	39
HDAC1 Inhibitor Screening Assay Kit	39
HDAC8 Inhibitor Screening Assay Kit	4
15(S)-HETE ELISA Kit	40
HIF-1α Transcription Factor Assay Kit	4
8-Hydroxyguanosine Monoclonal ELISA Kit	4
JC-1 Mitochondrial Membrane Potential Assay Kit	4
Lipid Hydroperoxide (LPO) Assay Kit	4
Lipid Hydroperoxide (LPO) Assay Kit (96 well)	4
LSD1 Inhibitor Screening Assay Kit	4
Malachite Green Phosphate Assay Kit	48
Methyltransferase Colorimetric Assay Kit	4
Methyltransferase Fluorometric Assay Kit	4
MLL1 SAM-Screener™ Assay Kit	
Multi-Parameter Apoptosis Assay Kit	
Nitrate/Nitrite Colorimetric Assay Kit	5
Oct4 Transcription Factor Assay Kit	
p53 Transcription Factor Assay Kit	5

PAD4 Inhibitor Screening Assay Kit (Ammonia)	56
cPLA <sub>2</sub> Assay Kit	57
PPARδ Transcription Factor Assay Kit	60
Prolactin (human) ELISA Kit	61
Protein Carbonyl Colorimetric Assay Kit	65
Protein Determination Kit	65
SET7/9 SAM-Screener™ Assay Kit	68
SIRT1 Direct Fluorescent Screening Assay Kit	68
SIRT1 FRET-Based Screening Assay Kit	68
SIRT2 Direct Fluorescent Screening Assay Kit	
SIRT3 Direct Fluorescent Screening Assay Kit	69
Sulfenylated Protein Cell-Based Detection Kit	
Superoxide Dismutase Assay Kit	70
TBARS Assay Kit	71
TBARS (TCA Method) Assay Kit	71



#### Cardiovascular Disease

Aconitase Assay Kit	15
Alanine Transaminase Colorimetric Activity Assay Kit	16
Angiotensin II EIA Kit	17
ApoA1 (human) ELISA Kit	18
Arginine Vasopressin ELISA Kit	18
Ascorbate Assay Kit	18
Atriopeptin (rat) EIA Kit	18
Cholesterol Fluorometric Assay Kit	23
COX Activity Assay Kit	24
COX Colorimetric Inhibitor Screening Assay Kit	24
COX (ovine) Colorimetric Inhibitor Screening Assay Kit	25
COX Fluorescent Activity Assay Kit	25
COX Fluorescent Inhibitor Screening Assay Kit	25
COX (ovine) Inhibitor Screening Assay Kit	25
COX (ovine/human) Inhibitor Screening Assay Kit	25
COX-1 (human) Inhibitor Screening Assay Kit	25
COX-1 (ovine) Inhibitor Screening Assay Kit	26
COX-2 (human) Inhibitor Screening Assay Kit	26
COX-2 (ovine) Inhibitor Screening Assay Kit	26
C-Reactive Protein (human) ELISA Kit	26
Creatinine (serum) Colorimetric Assay Kit	26
Creatinine (urinary) Colorimetric Assay Kit	26
Endothelin ELISA Kit	30
Hemoglobin Colorimetric Assay Kit	40
HIF-1α Transcription Factor Assay Kit	40
LDL Uptake Cell-Based Assay Kit	45
Lipase Activity Assay Kit	47
LXR (human) Reporter Assays Panel	48
LXRα (human) Reporter Assay Kit	48
LXRβ (human) Reporter Assay Kit	48
Malachite Green Phosphate Assay Kit	48
Myeloperoxidase Inhibitor Screening Assay Kit	52
Myeloperoxidase Peroxidation Fluorometric Assay Kit	52
PAF Acetylhydrolase Assay Kit	56
PAF Acetylhydrolase Inhibitor Screening Assay Kit	57
6-keto Prostaglandin F <sub>1α</sub> ELISA Kit	63
Prostaglandin I Metabolite ELISA Kit	64
Protein Carbonyl Colorimetric Assay Kit	65

Protein Determination Kit	.65
Renin Inhibitor Screening Assay Kit	
Soluble Epoxide Hydrolase Inhibitor Screening Assay Kit	.69
SREBP-2 Transcription Factor Assay Kit	.70
Thromboxane B <sub>2</sub> ELISA Kit	.72
Thromboxane B <sub>2</sub> Express ELISA Kit - Monoclonal	.73
11-dehydro Thromboxane B. FLISA Kit - Monoclonal	73



### Cell Biology

7-AAD/CFSE Cell-Mediated Cytotoxicity Assay Kit	. 1
AcSDKP EIA Kit	. 1
Adipogenesis Assay Kit	. 1
Adipolysis Assay Kit	1
Annexin V FITC Assay Kit	1
Autophagy/Cytotoxicity Dual Staining Kit	1
Autophlux Kit	1
Calcium Assay Kit	2
Caspase-3 Fluorescence Assay Kit	2
Cell Cycle Phase Determination Kit	2
CFSE Cell Division Assay Kit	2
CREB (Phospho-Ser <sup>133</sup> ) Transcription Factor Assay Kit	2
Cyclic AMP ELISA Kit	2
Cyclic AMP ELISA Kit (without Acetic Anhydride)	2
Cyclic AMP Select ELISA Kit	2
Cyclic GMP ELISA Kit	2
Cyclic GMP ELISA Kit (without Acetic Anhydride)	2
DNA Laddering Kit	7
DUB Activity Assay Kit	3
S-Glutathionylated Protein Detection Kit	3
Goat Anti-Human Fluorophore Selection Kit	3
Goat Anti-Mouse Fluorophore Selection Kit	3
Goat Anti-Rabbit Fluorophore Selection Kit	3
Hemoglobin Colorimetric Assay Kit	.4
HIF-1α Transcription Factor Assay Kit	.4
His-Tag Detection ELISA Kit	
JC-1 Mitochondrial Membrane Potential Assay Kit	.4
LC3 Interact Kit	.4
LDH Cytotoxicity Assay Kit	.4
LDL Uptake Cell-Based Assay Kit	.4
Live Cell Multiplex Assay Kit	.4
Malachite Green Phosphate Assay Kit	.4
MitoCheck® Citrate Synthase Activity Assay Kit	.4
MitoCheck® Complex I Activity Assay Kit	.4
MitoCheck® Complex II Activity Assay Kit	.4
MitoCheck® Complex II/III Activity Assay Kit	5
MitoCheck® Complex IV Activity Assay Kit	
MitoCheck® Complex V Activity Assay Kit	.5
MitoCheck® Mitochondrial (Tissue) Isolation Kit	7
MitoCheck® Mitochondrial OCR Assay Kit	.5
MTT Cell Proliferation Assay Kit	
Multi-Drug Resistance Assay Kit (Calcein AM)	5
Multi-Parameter Apoptosis Assay Kit	.5
Nuclear Extraction Kit	
Oct4 Transcription Factor Assay Kit	
Oxygen Consumption/Glycolysis Dual Assay Kit	5

Dxygen Consumption/MitoMembrane Potential Dual Assay Kit	55
PLA <sub>2</sub> (human Type IIA) ELISA Kit	58
OS Proteasome Assay Kit	64
Prostaglandin Screening ELISA Kit	64
Protein Determination Kit	65
Protein Synthesis Assay Kit	65
Pyruvate Assay Kit	65
Rabbit Anti-Goat Fluorophore Selection Kit	65
RARα (human) Reporter Assay Kit	65
RARβ (human) Reporter Assay Kit	66
RARγ (human) Reporter Assay Kit	66
RORγ (human) Reporter Assay Kit	67
RXR (human) Reporter Assays Panel	67
RXRα (human) Reporter Assay Kit	67
RXRβ (human) Reporter Assay Kit	67
RXRγ (human) Reporter Assay Kit	67
Secreted Alkaline Phosphatase Reporter Gene	
Assay Kit (Luminescence)	68
Sulfenylated Protein Cell-Based Detection Kit	70
SureLight® APC Conjugation Kit 1 mg IgG	71
SureLight® APC Conjugation Kit 2 mg IgG	71
SureLight® R-PE Conjugation Kit	71
MRE Mitochondrial Membrane Potential Assay Kit	73
Jbiquitin Interact Kit	74
Jbiquitin Link Kit	74
Jbiquitinated Protein Capture Kit	74
VST-1 Cell Proliferation Assay Kit	76
VST-8 Cell Proliferation Assay Kit	76
ATT Call Dealiferation Association	7/



#### Eicosanoids & Lipids

DX Activity Assay Kit	24
OX Colorimetric Inhibitor Screening Assay Kit	24
OX (ovine) Colorimetric Inhibitor Screening Assay Kit	25
DX Fluorescent Activity Assay Kit	25
OX Fluorescent Inhibitor Screening Assay Kit	25
OX (ovine) Inhibitor Screening Assay Kit	25
OX (ovine/human) Inhibitor Screening Assay Kit	25
DX-1 (human) Inhibitor Screening Assay Kit	25
DX-1 (ovine) Inhibitor Screening Assay Kit	26
DX-2 (human) Inhibitor Screening Assay Kit	26
DX-2 (ovine) Inhibitor Screening Assay Kit	26
steinyl Leukotriene ELISA Kit	28
steinyl Leukotriene Express ELISA Kit	
Receptor (rat) Reporter Assay Kit	30
<sup>7</sup> <sub>4</sub> Receptor (rat) Activation Assay Kit (cAMP)	30
AH Inhibitor Screening Assay Kit	32
BP4 Inhibitor/Ligand Screening Assay Kit	
ıprostenol ELISA Kit	33
ee Fatty Acid Fluorometric Assay Kit	
(S)-HETE ELISA Kit	40
Isoprostane Affinity Purification Kit	
Isoprostane ELISA Kit	
Isoprostane Express ELISA Kit	44

continued on next page

Latanoprost ELISA Kit	45
LDL Uptake Cell-Based Assay Kit	45
Leukotriene B <sub>4</sub> ELISA Kit	46
Leukotriene B <sub>4</sub> Express ELISA Kit	46
Leukotriene C <sub>4</sub> ELISA Kit	46
Leukotriene E <sub>4</sub> ELISA Kit	
Lipid Hydroperoxide (LPO) Assay Kit	47
Lipid Hydroperoxide (LPO) Assay Kit (96 well)	
Lipoxygenase Inhibitor Screening Assay Kit	47
Malachite Green Phosphate Assay Kit	
PAF Acetylhydrolase Assay Kit	
PAF Acetylhydrolase Inhibitor Screening Assay Kit	
tetranor-PGDM ELISA Kit	
Phosphatidylcholine Colorimetric Assay Kit	
cPLA <sub>2</sub> Assay Kit	
sPLA <sub>2</sub> Assay Kit	
sPLA <sub>2</sub> (human Type IIA) ELISA Kit	
sPLA <sub>2</sub> (Type V) Inhibitor Screening Assay Kit	
Prostaglandin D Synthase (hematopoietic-type) FP-Based	
Inhibitor Screening Assay Kit - Green	
Prostaglandin D <sub>2</sub> ELISA Kit	
Prostaglandin D <sub>2</sub> Express ELISA Kit	
Prostaglandin D <sub>2</sub> -MOX ELISA Kit	
Prostaglandin D <sub>2</sub> -MOX Express ELISA Kit	
Prostaglandin E Metabolite ELISA Kit	
Prostaglandin E <sub>2</sub> ELISA Kit - Monoclonal	
Prostaglandin E <sub>2</sub> Express ELISA Kit	
6-keto Prostaglandin F <sub>1α</sub> ELISA Kit	
Prostaglandin F <sub>2α</sub> ELISA Kit	
11β-Prostaglandin F <sub>2α</sub> ELISA Kit	
13,14-dihydro-15-keto Prostaglandin F <sub>2a</sub> ELISA Kit	
Prostaglandin I Metabolite ELISA Kit	
Prostaglandin Screening ELISA Kit	
Protein Determination Kit	
Resolvin D1 ELISA Kit	
Resolvin D2 ELISA Kit	
Soluble Epoxide Hydrolase Cell-Based Assay Kit	
Sphingomyelin Colorimetric Assay Kit	
STAT-8-Isoprostane ELISA Kit	
Thromboxane B <sub>2</sub> ELISA Kit	
Thromboxane B <sub>2</sub> Express ELISA Kit - Monoclonal	
11-dehydro Thromboxane B <sub>2</sub> ELISA Kit - Monoclonal	73

# Epigenetics

BAZ2B bromodomain TR-FRET Assay Kit	19
RD2 bromodomain 1 TR-FRET Assay Kit	19
RD2 bromodomain 2 TR-FRET Assay Kit	19
RD2 bromodomains 1 and 2 TR-FRET Assay Kit	20
RD3 bromodomain 1 TR-FRET Assay Kit	20
RD3 bromodomain 2 TR-FRET Assay Kit	20
RD3 bromodomains 1 and 2 TR-FRET Assay Kit	20
RD4 bromodomain 1 TR-FRET Assay Kit	20
RD4 bromodomain 2 TR-FRET Assay Kit	20
RD4 bromodomains 1 and 2 TR-FRET Assay Kit	21
RDT bromodomain 1 TR-FRET Assay Kit	21

BRG1 bromodomain TR-FRET Assay Kit21	L
BRM bromodomain TR-FRET Assay Kit21	L
CBP bromodomain TR-FRET Assay Kit22	)
DNA Methylation ELISA Kit29	)
EpiTitan™ Histone Peptide Array31	L
GLP SAM-Screener™ Assay Kit35	5
HAT Inhibitor Screening Assay Kit39	)
HDAC Cell-Based Activity Assay Kit39	)
HDAC Fluorometric Activity Assay Kit39	)
HDAC1 Inhibitor Screening Assay Kit39	)
HDAC8 Inhibitor Screening Assay Kit40	)
JMJD2A Tudor Domains TR-FRET Assay Kit44	1
L3MBTL1 MBT Domains TR-FRET Assay Kit44	1
LSD1 Inhibitor Screening Assay Kit47	7
Methyltransferase Colorimetric Assay Kit49	)
Methyltransferase Fluorometric Assay Kit49	)
MLL1 SAM-Screener™ Assay Kit50	)
SET7/9 SAM-Screener™ Assay Kit68	3
SIRT1 Direct Fluorescent Screening Assay Kit68	3
SIRT1 FRET-Based Screening Assay Kit68	3
SIRT2 Direct Fluorescent Screening Assay Kit68	3
SIRT3 Direct Fluorescent Screening Assay Kit69	)
TAF1 bromodomain 1 TR-FRET Assay Kit71	L
TAF1 bromodomains 1 and 2 TR-FRET Assay Kit71	L

# Immunology & Inflammation

7-AAD/CFSE Cell-Mediated Cytotoxicity Assay Kit	15
Anti-Citrullinated Human Fibrinogen Assay Kit (mouse)	17
Anti-Ovalbumin IgE (mouse) ELISA Kit	17
Anti-Ovalbumin IgG, (mouse) ELISA Kit	17
Autotaxin Inhibitor Screening Assay Kit	19
CFSE Cell Division Assay Kit	23
COX Activity Assay Kit	24
COX Colorimetric Inhibitor Screening Assay Kit	24
COX (ovine) Colorimetric Inhibitor Screening Assay Kit	25
COX Fluorescent Activity Assay Kit	25
COX Fluorescent Inhibitor Screening Assay Kit	25
COX (ovine) Inhibitor Screening Assay Kit	25
COX (ovine/human) Inhibitor Screening Assay Kit	25
COX-1 (human) Inhibitor Screening Assay Kit	25
COX-1 (ovine) Inhibitor Screening Assay Kit	26
COX-2 (human) Inhibitor Screening Assay Kit	26
COX-2 (ovine) Inhibitor Screening Assay Kit	26
C-Reactive Protein (human) ELISA Kit	26
Cysteinyl Leukotriene ELISA Kit	28
Cysteinyl Leukotriene Express ELISA Kit	28
EP <sub>2</sub> Receptor (rat) Reporter Assay Kit	30
EP <sub>4</sub> Receptor (rat) Activation Assay Kit (cAMP)	30
EP <sub>4</sub> Receptor (rat) Reporter Assay Kit	31
15(S)-HETE ELISA Kit	40
Histamine EIA Kit	40
Human Therapeutic IgG1 ELISA Kit	41
Human Therapeutic IgG2 ELISA Kit	
Human Therapeutic IgG4 ELISA Kit	41
Hydrogen Peroxide Cell-Based Assay Kit	41

Interleukin-1α (human) ELISA Kit	43
Interleukin-1β (human) ELISA Kit	43
Interleukin-6 (human) ELISA Kit	43
Interleukin-6 (mouse) ELISA Kit	43
LDH Cytotoxicity Assay Kit	45
Leukotriene B <sub>4</sub> ELISA Kit	46
Leukotriene B <sub>4</sub> Express ELISA Kit	46
Leukotriene C <sub>4</sub> ELISA Kit	46
Leukotriene E <sub>4</sub> ELISA Kit	46
Malachite Green Phosphate Assay Kit	48
Mouse Anti-Type II Collagen IgG Assay Kit (bovine)	51
Mouse IgG ELISA Kit	51
Myeloperoxidase Chlorination Fluorometric Assay Kit	52
Myeloperoxidase Inhibitor Screening Assay Kit	52
Myeloperoxidase Peroxidation Fluorometric Assay Kit	52
Neutrophil Elastase Activity Assay Kit	52
Neutrophil Extracellular Trap (MPO-Histone) ELISA Kit	52
Neutrophil Extracellular Trap (NET) Assay Kit	53
Neutrophil (mouse) Isolation Kit	77
Neutrophil/Monocyte Respiratory Burst Assay Kit	53
Neutrophil Myeloperoxidase Activity Assay Kit	53
NF-ĸB (human p50) Transcription Factor Assay Kit	53
NF-ĸB (p65) Transcription Factor Assay Kit	53
Nrf2 (human) Reporter Assay Kit	54
PAD4 Autoantibody ELISA Kit	56
PAD4 Inhibitor Screening Assay Kit (Ammonia)	56
PAF Acetylhydrolase Inhibitor Screening Assay Kit	57
tetranor-PGDM ELISA Kit	57
Phagocytosis Assay Kit (IgG FITC)	57
Phagocytosis Assay Kit (IgG PE)	
cPLA <sub>2</sub> Assay Kit	57
sPLA <sub>2</sub> (human Type IIA) ELISA Kit	58
PPARδ Transcription Factor Assay Kit	60
Prostaglandin D Synthase (hematopoietic-type) FP-Based	
Inhibitor Screening Assay Kit - Green	
Prostaglandin D <sub>2</sub> ELISA Kit	
Prostaglandin D <sub>2</sub> Express ELISA Kit	
Prostaglandin D <sub>2</sub> -MOX ELISA Kit	62
Prostaglandin E Metabolite ELISA Kit	
Prostaglandin E <sub>2</sub> ELISA Kit - Monoclonal	
Prostaglandin E <sub>2</sub> Express ELISA Kit	63
11β-Prostaglandin $F_{2\alpha}$ ELISA Kit	63
Prostaglandin Screening ELISA Kit	
Protein Carbonyl Colorimetric Assay Kit	65
Protein Determination Kit	
Resolvin D1 ELISA Kit	66
Resolvin D2 ELISA Kit	
Soluble Epoxide Hydrolase Inhibitor Screening Assay Kit	
Thromboxane B <sub>2</sub> ELISA Kit	
Thromboxane B <sub>2</sub> Express ELISA Kit - Monoclonal	
11-dehydro Thromboxane B <sub>2</sub> ELISA Kit - Monoclonal	
TNF-α (human) ELISA Kit	73

TNF-α (mouse) ELISA Kit.....



- A A	0+0	0	icia
IVI	eta	arai	usrr
	CLUI	$\sigma \sigma$	1311

cid Phosphatase Colorimetric Assay Kit	. 1
conitase Assay Kit	. 1
dipogenesis Assay Kit	. 1
dipolysis Assay Kit	. 1
diponectin (human) EIA Kit	. 1
diponectin (human) EIA Kit (HS)	. 1
diponectin (mouse) EIA Kit	
hR (human) Assay Kit	. 1
lanine Transaminase Colorimetric Activity Assay Kit	
ldehyde Dehydrogenase Activity Assay Kit	
ldosterone ELISA Kit	. 1
ntioxidant Assay Kit	. 1
poA1 (human) ELISA Kit	
IR (human) Reporter Assay Kit	
rginine Vasopressin ELISA Kit	. 1
scorbate Assay Kit	. 1
Calcium Assay Kit	.2
Cholesterol Cell-Based Detection Assay Kit	.2
Cholesterol Fluorometric Assay Kit	.2
Cholesterol Uptake Cell-Based Assay Kit	.2
Corticosterone ELISA Kit	.2
Cortisol ELISA Kit	.2
Cortisol Express ELISA Kit	
Creatinine (serum) Colorimetric Assay Kit	
reatinine (urinary) Colorimetric Assay Kit	
OPP (IV) Inhibitor Screening Assay Kit	
ndothelin ELISA Kit	.3
nterolactone ELISA Kit	
P <sub>4</sub> Receptor (rat) Reporter Assay Kit	
R (human) Reporter Assays Panel	
Rα (human) Reporter Assay Kit	
Rβ (human) Reporter Assay Kit	
RRα (human) Reporter Assay Kit	
RRγ (human) Reporter Assay Kit	
stradiol ELISA Kit	
striol ELISA Kit	
ABP4 (human) EIA Kit	
ABP4 Inhibitor/Ligand Screening Assay Kit	
luprostenol ELISA Kit	
ollicle-Stimulating Hormone ELISA Kit	
ree Fatty Acid Fluorometric Assay Kit	
XR (human) Reporter Assay Kit	
cylated Ghrelin (dog) Easy Sampling EIA Kit	
cylated Ghrelin (human) Easy Sampling EIA Kit	
cylated Ghrelin (human) Express EIA Kit	
cylated Ghrelin (mouse, rat) Easy Sampling EIA Kit	
cylated Ghrelin (mouse, rat) Express EIA Kit	
cylated Ghrelin (pig) EIA Kit	
Jnacylated Ghrelin (dog) Easy Sampling EIA Kit	
Jnacylated Ghrelin (human) Easy Sampling EIA Kit	
Jnacylated Ghrelin (mouse, rat) Easy Sampling EIA Kit	
Jnacylated Ghrelin (pig) EIA Kit	
Jnacylated Ghrelin (human) Express EIA Kit	
Inacylated Ghrelin (mouse, rat) Express EIA Kit	
continued on next n	no

Glucose Colorimetric Assay Kit	36
Glucose-6-Phosphate Dehydrogenase Activity Assay Kit	36
Glucose-6-Phosphate Fluorometric Assay Kit	36
Glucose Uptake Cell-Based Assay Kit	36
Glycerol Cell-Based Assay Kit	37
Glycerol Colorimetric Assay Kit	38
Glycogen Assay Kit	38
Glycolysis Cell-Based Assay Kit	38
GR (human) Reporter Assay Kit	39
Growth Hormone (rat) EIA Kit	39
β-Hydroxybutyrate (Ketone Body) Colorimetric Assay Kit	42
β-Hydroxybutyrate (Ketone Body) Fluorometric Assay Kit	42
20-Hydroxyecdysone EIA Kit	42
Insulin (rat) EIA Kit	42
D-Lactate Assay Kit	44
L-Lactate Assay Kit	45
Leptin (human) EIA Kit	45
Leptin (mouse/rat) EIA Kit	46
Leptin Receptor (human) EIA Kit	46
Lipase Activity Assay Kit	47
Lipid Droplets Fluorescence Assay Kit	47
Lipid Hydroperoxide (LPO) Assay Kit	47
Lipid Hydroperoxide (LPO) Assay Kit (96 well)	47
Lipoxygenase Inhibitor Screening Assay Kit	47
Luteinizing Hormone ELISA Kit	48
Malachite Green Phosphate Assay Kit	48
Melanocortin-3 Receptor Reporter Assay Kit	48
Melanocortin-4 Receptor Reporter Assay Kit	48
2-Methoxyestradiol ELISA Kit	49
MitoCheck® Citrate Synthase Activity Assay Kit	49
MitoCheck® Complex I Activity Assay Kit	49
MitoCheck® Complex II Activity Assay Kit	49
MitoCheck® Complex II/III Activity Assay Kit	50
MitoCheck® Complex IV Activity Assay Kit	
MitoCheck® Complex V Activity Assay Kit	50
MitoCheck® Mitochondrial (Tissue) Isolation Kit	77
MitoCheck® Mitochondrial OCR Assay Kit	50
MR (human) Reporter Assay Kit	51
NAD/NADH Cell-Based Assay Kit	52
Orexin 1 Receptor Reporter Assay Kit	
Oxygen Consumption Rate Assay Kit (MitoXpress $^{\mathbb{R}}\text{-Xtra HS Method})$	55
Oxygen Consumption/Glycolysis Dual Assay Kit	55
Oxygen Consumption/MitoMembrane Potential Dual Assay Kit	55
Oxytocin ELISA Kit	56
PAF Acetylhydrolase Assay Kit	56
PAF Acetylhydrolase Inhibitor Screening Assay Kit	57
PGR (human) Reporter Assay Kit	57
Phosphatidylcholine Colorimetric Assay Kit	57
PPAR (human) Reporter Assays Panel	58
PPARα (human) Reporter Assay Kit	58
PPARα (mouse) Reporter Assay Kit	58
PPARα (rat) Reporter Assay Kit	
PPARα Transcription Factor Assay Kit	59
PPAR $\alpha$ , $\delta$ , $\gamma$ Complete Transcription Factor Assay Kit	
PPARδ (human) Reporter Assay Kit	
PPARδ (mouse) Reporter Assay Kit	59
PPARδ (rat) Reporter Assay Kit	
PPARδ Transcription Factor Assay Kit	60

PPARγ (human) Reporter Assay Kit	60
PPARγ (mouse/rat) Reporter Assay Kit	60
PPARγ Transcription Factor Assay Kit	60
Progesterone ELISA Kit	60
Prolactin (human) ELISA Kit	61
Prolactin (rat) EIA Kit	61
Prostaglandin E Metabolite ELISA Kit	62
Prostaglandin E <sub>2</sub> ELISA Kit - Monoclonal	62
Prostaglandin E <sub>2</sub> Express ELISA Kit	63
Prostaglandin F <sub>2α</sub> ELISA Kit	63
L3,14-dihydro-15-keto Prostaglandin F <sub>2a</sub> ELISA Kit	
L7-phenyl trinor Prostaglandin F <sub>2a</sub> ELISA Kit	
Protein Determination Kit	
Pyruvate Assay Kit	65
RAR (human) Reporter Assays Panel	
Renin Inhibitor Screening Assay Kit	66
Resistin (human) EIA Kit	
Resistin (rat) EIA Kit	
SIRT1 Direct Fluorescent Screening Assay Kit	
SIRT1 FRET-Based Screening Assay Kit	
SIRT2 Direct Fluorescent Screening Assay Kit	68
SIRT3 Direct Fluorescent Screening Assay Kit	
Soluble Epoxide Hydrolase Inhibitor Screening Assay Kit	
SREBP-1 Transcription Factor Assay Kit	
SREBP-2 Cell-Based Translocation Assay Kit	69
SREBP-2 Transcription Factor Assay Kit	
Steatosis Colorimetric Assay Kit	70
Substance P ELISA Kit	7C
FBARS Assay Kit	71
FBARS (TCA Method) Assay Kit	71
Testosterone ELISA Kit	71
l 1-keto Testosterone ELISA Kit	72
Thyroid-Stimulating Hormone ELISA Kit	73
TMRE Mitochondrial Membrane Potential Assay Kit	73
FR (human) Reporter Assays Panel	74
ΓRα (human) Reporter Assay Kit	
ΓRβ (human) Reporter Assay Kit	74
Friglyceride Colorimetric Assay Kit	74
Jrea Fluorometric Assay Kit	74
Jric Acid Assay Kit	
/DR (human) Reporter Assay Kit	
/itamin D ELISA Kit	

# > Neuroscience

CGRP (human) EIA Kit	23
CGRP (rat) EIA Kit	23
EP <sub>4</sub> Receptor (rat) Activation Assay Kit (cAMP)	30
FAAH Inhibitor Screening Assay Kit	32
Fluprostenol ELISA Kit	33
GFAP (human) EIA Kit	33
Latanoprost ELISA Kit	45
Malachite Green Phosphate Assay Kit	48
Melanocortin-3 Receptor Reporter Assay Kit	48
Melanocortin-4 Receptor Reporter Assay Kit	48
Monoacylglycerol Lipase Inhibitor Screening Assay Kit	50

Orexin 1 Receptor Reporter Assay Kit	55
Orexin 2 Receptor Reporter Assay Kit	55
tetranor-PGDM ELISA Kit	57
Prostaglandin D <sub>2</sub> ELISA Kit	61
Prostaglandin D <sub>2</sub> Express ELISA Kit	61
Prostaglandin D <sub>2</sub> -MOX ELISA Kit	62
Prostaglandin D <sub>2</sub> -MOX Express ELISA Kit	62
11β-Prostaglandin F <sub>2α</sub> ELISA Kit	63
Protein Determination Kit	
Resolvin D1 ELISA Kit	66
Substance P ELISA Kit	70

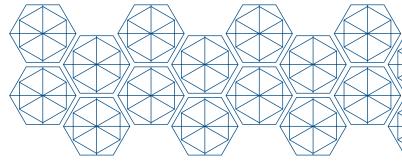


# Nitric Oxide & Oxidative Injury

Antioxidant Assay Kit	1
Catalase Assay Kit	2
Catalase Assay Kit (without Hydrogen Peroxide)	2
DHN-MA EIA Kit	2
DNA/RNA Oxidative Damage ELISA Kit	2
Glutaredoxin Fluorescent Activity Assay Kit	3
Glutathione Assay Kit	3
Glutathione Cell-Based Detection Kit (Blue Fluorescence)	3
Glutathione Peroxidase Assay Kit	3
Glutathione Reductase Assay Kit	3
Glutathione S-Transferase Assay Kit	3
S-Glutathionylated Protein Detection Kit	3
Glyco-SPOT DNA Repair Assay Kit	3
8-Hydroxyguanosine Monoclonal ELISA Kit	4
8-Isoprostane Affinity Purification Kit	7
8-Isoprostane ELISA Kit	4
8-Isoprostane Express ELISA Kit	4
Lipid Hydroperoxide (LPO) Assay Kit	4
Lipid Hydroperoxide (LPO) Assay Kit (96 well)	4
Malachite Green Phosphate Assay Kit	4
Nitrate/Nitrite Colorimetric Assay Kit	5
Nitrate/Nitrite Colorimetric Assay Kit (LDH method)	5
Nitrate/Nitrite Fluorometric Assay Kit	5
S-Nitrosylated Protein Detection Kit	5
NOS Activity Assay Kit	5
Nrf2 Transcription Factor Assay Kit	5
Protein Carbonyl Colorimetric Assay Kit	6
Protein Determination Kit	6
STAT-8-Isoprostane ELISA Kit	7
Sulfenylated Protein Cell-Based Detection Kit	
Superoxide Dismutase Assay Kit	7
TBARS Assay Kit	7
TBARS (TCA Method) Assay Kit	7
Thiol Detection Assay Kit	7
Thioredoxin Activity Fluorescent V2 Assay Kit	
Thioredoxin Reductase Colorimetric Assay Kit	7
Thioredoxin Reductase Fluorescent Activity Assay Kit	7
Thioredoxin/Thioredoxin Reductase Mammalian Assay Kit	
Xanthine Oxidase Fluorometric Assay Kit	7



Manine transaminase Coloninetric Activity Assay Kit	TC
Aldehyde Dehydrogenase Activity Assay Kit	16
ASP ELISA Kit	18
CAR1 (human) Reporter Assay Kit	21
CAR3 (human) Reporter Assay Kit	22
CYP1A1/2 Induction Reporter Assay Kit	27
CYP2B6 Induction Reporter Assay Kit	28
CYP2C9 Induction Reporter Assay Kit	28
Human Serum Albumin ELISA Kit	41
Human Therapeutic IgG1 ELISA Kit	41
Human Therapeutic IgG2 ELISA Kit	41
Human Therapeutic IgG4 ELISA Kit	41
DH Cytotoxicity Assay Kit	45
Malachite Green Phosphate Assay Kit	48
MitoCheck® Citrate Synthase Activity Assay Kit	49
الله المالة المالة AitoCheck® Complex ا Activity Assay Kit	49
AitoCheck® Complex II Activity Assay Kit	49
NitoCheck® Complex II/III Activity Assay Kit	50
AitoCheck® Complex IV Activity Assay Kit	50
MitoCheck® Complex V Activity Assay Kit	50
AitoCheck® Mitochondrial (Tissue) Isolation Kit	77
AitoCheck® Mitochondrial OCR Assay Kit	50
MTT Cell Proliferation Assay Kit	51
Multi-Drug Resistance Assay Kit (Calcein AM)	51
Oxygen Consumption Rate Assay Kit (MitoXpress®-Xtra HS Method)	55
Oxygen Consumption/Glycolysis Dual Assay Kit	55
Oxygen Consumption/MitoMembrane Potential Dual Assay Kit	
53 Transcription Factor Assay Kit	
Pregnane X (human) Receptor Assay Kit	60
Pregnane X (rat) Receptor Assay Kit	
Protein Determination Kit	65
Semi-Quantitative Biomarker ELISA Component Kit (anti-mouse)	68
Semi-Quantitative Biomarker ELISA Component Kit (anti-rabbit)	68
steatosis Colorimetric Assay Kit	70
/itellogenin (carp) ELISA Kit	75
/itellogenin (fathead minnow) ELISA Kit	
/itellogenin (medaka) ELISA Kit	75
/itellogenin (rainbow trout) ELISA Kit	75
/itellogenin (salmonid) Semi-Quantitative ELISA Kit	
/itellogenin (zebrafish) ELISA Kit	
VST-1 Cell Proliferation Assay Kit	
VST-8 Cell Proliferation Assay Kit	76
KTT Cell Proliferation Assay Kit	



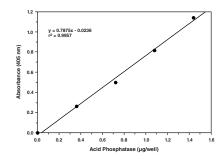
### KIT INFORMATION (ALPHABETICAL)

#### 7-AAD/CFSE Cell-Mediated Cytotoxicity Assay Kit 600120

- Label target cells with highly fluorescent 5-(6)-carboxyfluorescein diacetate succinimidyl ester (CFSE) for easy identification
- Dead cells stain with 7-amino actinomycin D (7-AAD)
- Analyze by flow cytometry to assess target cell killing

#### Acid Phosphatase Colorimetric Assay Kit 10008051

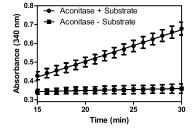
- Measure acid phosphatase (AP) activity in plasma, serum, urine, and
- Assay 44 samples in duplicate/96-well plate
- Includes L-tartrate for measuring tartrate resistant AP activity
- Plate-based colorimetric measurement (405-414 nm)



#### Aconitase Assay Kit

705502

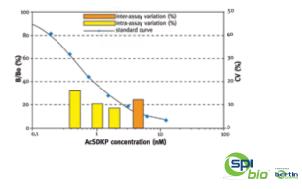
- Measure aconitase activity in tissue homogenates or cell lysates
- Assay 23 samples in duplicate
- Measure aconitase activity down to 1.7 U/ml
- Plate-based colorimetric measurement (340 nm)



#### AcSDKP EIA Kit

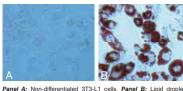
589451

- Measure N-Acetyl Ser-Asp-Lys-Pro (AcSDKP) in mammalian plasma and urine samples
- Assay 33 samples in duplicate
- Incubation: Overnight



#### Adipogenesis Assay Kit

- Contains the reagents needed for induction of adipogenesis in vitro
- Contains the reagents needed to stain lipid droplets in adipocytes
- Assess by microscopy or plate reader



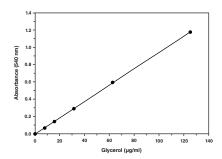
accumulate upon differentiation of 3T3-L1 cells, visualized by Oil Red O Solution staining.

#### Adipolysis Assay Kit

10009381

10006908

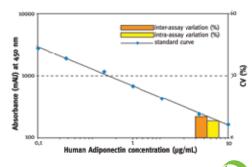
- Screen compounds involved in lipid storage and metabolism
- Contains the reagents needed for induction of adipogenesis in vitro, to release glycerol from stored lipid droplets, and to accurately measure the released glycerol



#### Adiponectin (human) EIA Kit

500641

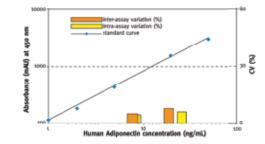
- Measure adiponectin in human serum and plasma samples
- Assay 40 samples in duplicate
- Incubation: 2 hours





#### Adiponectin (human) EIA Kit (HS)

- Measure adiponectin in human serum, plasma, breast milk, urine, and tissue culture supernatant samples
- Assay 41 samples in duplicate
- Incubation: 2 x 1 hour

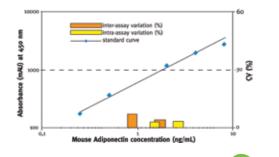


#### Adiponectin (mouse) EIA Kit

10007620

10007619

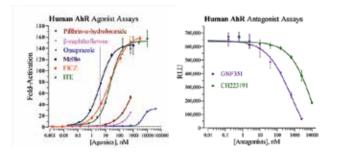
- Measure adiponectin in mouse samples
- Assay 41 samples in duplicate
- Incubations: 2 x 1 hour | 30 minutes



#### AhR (human) Assay Kit

16523

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

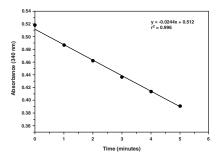




#### Alanine Transaminase Colorimetric **Activity Assay Kit**

700260

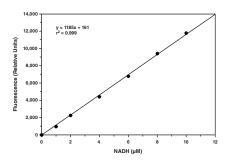
- Measure alanine transaminase (ALT) activity in serum, plasma, tissue samples, and cell lysates
  - Assay 47 samples in duplicate
  - Measure ALT activity down to 0.006 U/ml
  - Plate-based colorimetric measurement (340 nm)



#### Aldehyde Dehydrogenase Activity Assay Kit

700800

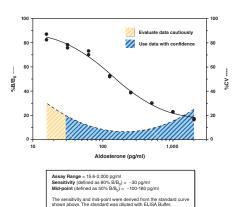
- Measure aldehyde dehydrogenase (ALDH) activity in tissue homogenates, cell culture samples, and purified ALDH preparations
- Assay 20 samples in duplicate
- Measure ALDH activity down to 0.4 U/ml
- Plate-based fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### Aldosterone ELISA Kit

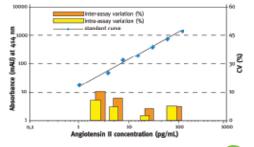
501090

- Measure aldosterone, the principle mineralocorticoid controlling sodium and potassium balance
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure aldosterone levels down to 30 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



WWW.**Caymanchem**.com Kit Information (Listed Alphabetically) **(**16)

- Measure angiontensin II in mammalian plasma, serum, and cell culture supernatant samples
- Assay 33 samples in duplicate
- Incubations: 1 hour | Overnight
- Angiotensin II Inhibitor Cocktail (Item No. 9000681) and phenyl cartridges (Item No. 400048) available for sample storage and clean up



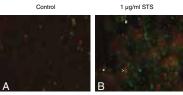


#### Annexin V FITC Assay Kit

600300

589301

- Identify cells undergoing apoptosis
- Assess by flow cytometry or fluorescence microscopy
- Sufficient reagent to test 500 samples
- 30 minute staining protocol

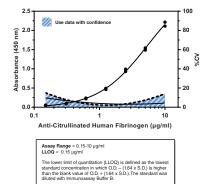


Staurosporine (STS) induced apoptosis in RAW 264.7 cells. Annexin V-FITC and Pl labeled cells were visualized with fluorescence microscopy. Annexin V-FITC positive cells (\*) are early apoptotic, double positive cells (>) are late apoptotic, and PI single positive cells

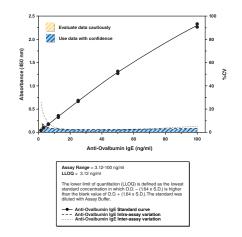
#### Anti-Citrullinated Human Fibrinogen Assay Kit (mouse)

501270

- Measure antibodies specific for citrullinated human fibrinogen
- Use with mouse plasma or serum
- Assay Range: 0.15-10 μg/ml with an LLOQ of 0.15 μg/ml
- Human fibrinogen affinity sorbent included for purification from any antibodies capable of reacting with non-citrullinated (unmodified) fibrinogen prior to analysis



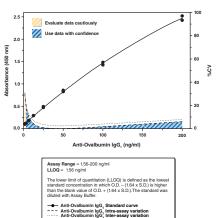
- Measure mouse anti-ovalbumin IgE (anti-OVA IgE) in mouse plasma, serum, or BAL fluids
- Employ a mouse anti-OVA IgE monoclonal antibody standard for reproducible results
- Assay Range: 3.12-100 ng/ml with an LLOQ of 3.12 ng/ml
- Rapid assay; get results in under 4 hours



#### Anti-Ovalbumin IgG<sub>1</sub> (mouse) ELISA Kit

500830

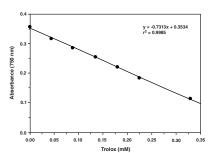
- Measure mouse anti-ovalbumin IgG, (anti-OVA IgG,) in mouse plasma, serum, or BAL fluids
- Employ an affinity-purified mouse anti-OVA IgG, polyclonal antibody standard for accurate results
- Assay Range: 1.56-200 ng/ml with an LLOQ of 1.56 ng/ml
- Rapid assay; get results in under 4 hours



#### Antioxidant Assay Kit

709001

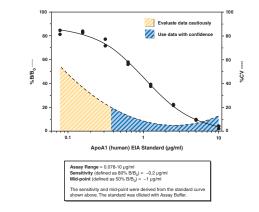
- Measure the total antioxidant capacity of plasma, serum, urine, saliva, or cell lysates
- Assay 41 samples in duplicate
- Measure antioxidant capacity in Trolox equivalents down to 44 μM
- Plate-based colorimetric measurement (750 or 405 nm)



#### • Measure human apolipoprotein A1 (ApoA1), a major component of HDL, in plasma or serum

- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure ApoA1 levels down to 200 ng/ml
- Rapid assay; get results in under 4 hours

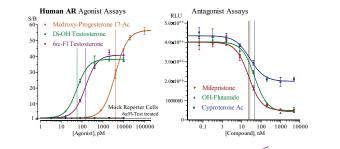
ApoA1 (human) ELISA Kit



#### AR (human) Reporter Assay Kit

15756

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

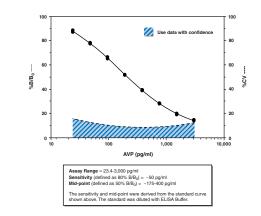


#### Arginine Vasopressin ELISA Kit

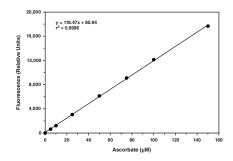
583951

**MINDIGO** 

- Measure arginine vasopressin (AVP), a nine amino acid peptide that regulates water secretion in all mammals
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure plasma AVP levels down to 50 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



- Quantify ascorbate from plasma, serum, urine, and fruit juices
- Assay 40 samples in duplicate
- Assay Range: 5-150 μM
- Plate-based fluorometric measurement (ex 340-350 nm, em 420-430 nm)



#### **ASP ELISA Kit**

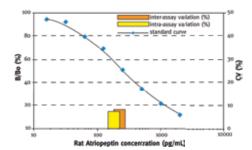
10008673

- A specific assay for domoic acid, a neurotoxin responsible for amnesic shellfish poisoning (ASP)
- See www.caymanchem.com for more details



#### Atriopeptin (rat) EIA Kit

- Measure atriopeptin in rat plasma or tissue extract samples
- Assay 33 samples in duplicate
- Incubations: 1 hour | Overnight
- Tissue extract should be purified using SPE catridges (C-18) (Item No. 400020)





Autophagy/Cytotoxicity Dual Staining Kit

• Propidium iodide included for dead cell exclusion

• Tamoxifen included as an autosome-inducing control

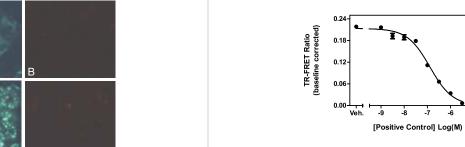
autophagic vacuoles

• Utilize the fluorescent reagent monodansylcadaverine to label

600140

BAZ2B bromodomain TR-FRET Assay Kit

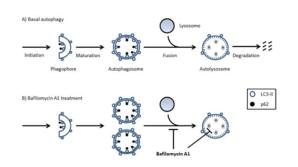
- 600710
- Homogeneous, time-resolved fluorescence energy transfer
- Plate-based, time-resolved fluorometric measurement (ex 340 nm,



#### Autophlux Kit

15980

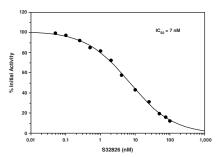
- Determine effect of your experiment on autophagy
- Identify and characterize autophagy modulators (activators/inhibitors)
- Assess result of gene knockdown (siRNA/shRNA) on autophagy
- Evaluate outcome of plasmid transfection/protein expression on
- Investigate autophagy in specific cell lines



#### Autotaxin Inhibitor Screening Assay Kit

700580

- Screen for inhibitors of autotaxin
- Assay 45 samples in duplicate
- Plate-based colorimetric measurment (405-415 nm)



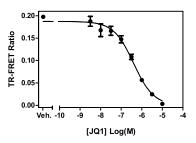
#### (TR-FRET) assay to screen for inhibitors • Designed for high-throughput screening

- em 620 and 670 nm)

#### BRD2 bromodomain 1 TR-FRET Assay Kit

600500

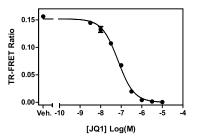
- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRD2 bromodomain 2 TR-FRET Assay Kit

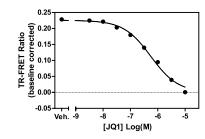
600510

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRD2 bromodomains 1 and 2 TR-FRET Assay Kit 600810

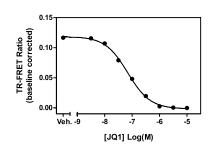
- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- · Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRD3 bromodomain 1 TR-FRET Assay Kit

600630

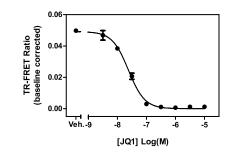
- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRD3 bromodomain 2 TR-FRET Assay Kit

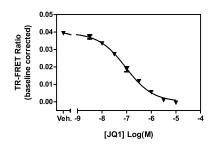
600640

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRD3 bromodomains 1 and 2 TR-FRET Assay Kit 600820

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)

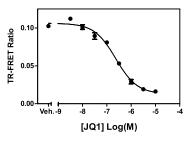


#### BRD4 bromodomain 1 TR-FRET Assay Kit

600520

600530

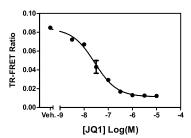
- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRD4 bromodomain 2 TR-FRET Assay Kit

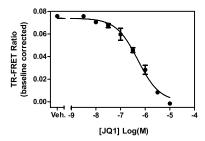
• Homogeneous, time-resolved fluorescence energy transfer

- (TR-FRET) assay to screen for inhibitors • Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



BRD4 bromodomains 1 and 2 TR-FRET Assay Kit 600830

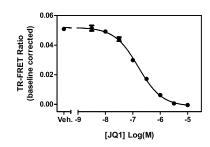
- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- · Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRDT bromodomain 1 TR-FRET Assay Kit

600650

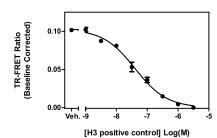
- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### BRG1 bromodomain TR-FRET Assay Kit

600720

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



BRM bromodomain TR-FRET Assay Kit

(TR-FRET) assay to screen for inhibitors

• Designed for high-throughput screening

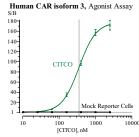
em 620 and 670 nm)

• Homogeneous, time-resolved fluorescence energy transfer

• Plate-based, time-resolved fluorometric measurement (ex 340 nm,

600730

- - See www.caymanchem.com for more details



# **MINDIGO**

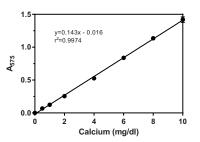
#### Calcium Assay Kit

701220

• Measure calcium in plasma, serum, urine, tissue homogenates, and cell lysates

[Positive Control] Log(M)

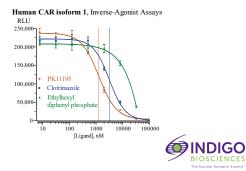
- Assay 40 samples in duplicate
- Measure calcium down to 0.25 mg/dl
- Plate-based, colorimetric measurement (570-590 nm)



#### CAR1 (human) Reporter Assay Kit

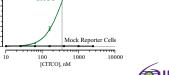
15748

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



#### CAR3 (human) Reporter Assay Kit

- A nuclear receptor cell-based reporter assay

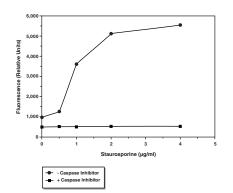


#### Caspase-3 Fluorescence Assay Kit

10009135

15747

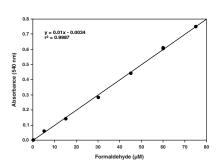
- Measure caspase-3 enzymatic activity levels in cell lysates
- Works with caspase-3 from multiple species
- Active caspase-3 generates a fluorescent product that is easily guantified



#### Catalase Assay Kit

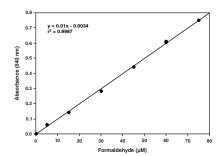
707002

- Measure catalase activity in plasma, serum, erythrocyte lysates, tissue homogenates, and cell lysates
- Assay 40 samples in duplicate
- Measure catalase activity down to 2 U/ml
- Plate-based, colorimetric measurement (540 nm)



#### Catalase Assay Kit (without Hydrogen Peroxide)

- Measure catalase activity in plasma, serum, erythrocyte lysates, tissue homogenates, and cell lysates
- Assay 40 samples in duplicate
- Measure catalase activity down to 2 U/ml
- Plate-based, colorimetric measurement (540 nm)

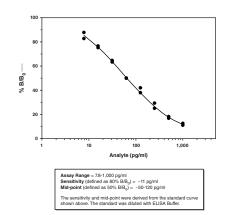


#### Cayman Practice ELISA Kit

10009658

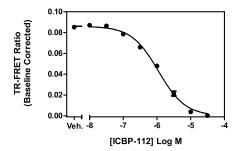
700910

- A practice ELISA kit for those unfamiliar with performing competitive ELISAs
- Easy to follow instructions with helpful illustrations
- Low cost



#### CBP bromodomain TR-FRET Assay Kit

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



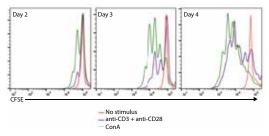
- Flow cytometric analysis of cell cycle progression (G<sub>0</sub>/G<sub>1</sub>, G<sub>2</sub>, or S phase)
- Contains reagents for fixing, permeabilizing, and staining cells
- Easy to follow protocol

#### **CFSE Cell Division Assay Kit**

10009853

10009349

- Flow cytometric analysis of cell proliferation
- Cell division results in a halving of the fluorescent signal; up to eight divisions can be monitored
- Compatible with staining for cell-surface and intracellular markers

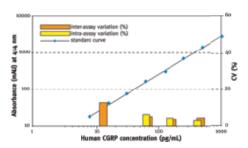


CD4+ T cells proliferate in response to T cell receptor antibodies and mitogenic stimuli

#### CGRP (human) EIA Kit

589101

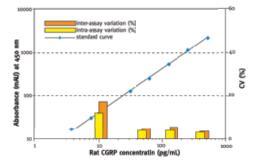
- Measure calcitonin gene-related peptide (CGRP) in human plasma, serum, CSF, bronchoalveolar lavage, cell extract, and cell culture supernatant samples
- Assay 33 samples in duplicate
- Incubation: Overnight
- SPE Catridges (C-18) (Item No. 400020) available for easy and safe extraction of CGRP from plasma samples





• Measure calcitonin gene-related peptide (CGRP) in mouse or rat plasma, serum, CSF, bronchoalveolar lavage, cell extract, and cell culture supernatant samples

- Assav 33 samples in duplicate
- Incubation: Overnight
- SPE Catridges (C-18) (Item No. 400020) available for easy and safe extraction of CGRP from plasma samples

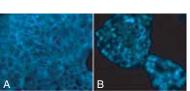




#### Cholesterol Cell-Based Detection Assay Kit

10009779

- Detect cholesterol uptake and localization within cultured cells
- Employ filipin III, a probe that fluoresces in the presence of cholesterol
- An inhibitor of cholesterol trafficking is included as a positive control

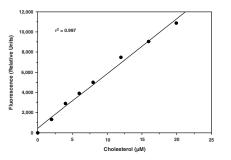


Accumulation of cholesterol in HepG2 cells in response to J-18666A for 48 hours. Panel A: Vehicle treated cells den

#### Cholesterol Fluorometric Assay Kit

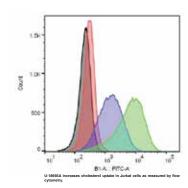
10007640

- Quantitation of total cholesterol in plasma or serum
- Assay 40 samples in duplicate
- Assay Range: 2-20 μM
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### • Detect cholesterol uptake within cultured cells by flow cytometry or plate reader

- Employ NBD cholesterol, a fluorescently-tagged cholesterol
- An inhibitor of cholesterol trafficking is included as a positive control
- The kit provides enough NBD cholesterol to test 250 samples

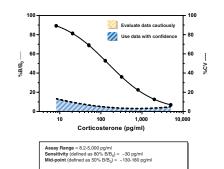


#### Corticosterone ELISA Kit

501320

600440

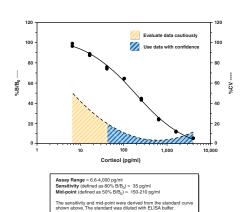
- Measure corticosterone in serum and rodent feces
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure corticosterone levels down to 30 pg/ml
- Incubation: Overnight | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



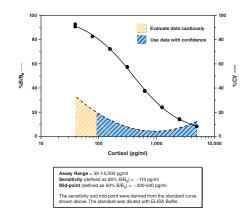
#### Cortisol ELISA Kit

500360

- Measure cortisol in urine, plasma, and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure cortisol levels down to 35 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



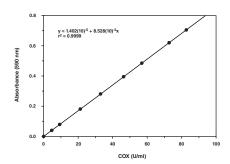
- Measure cortisol in urine, plasma, and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure cortisol levels down to 110 pg/ml
- Rapid assay; get results in under 4 hours



#### COX Activity Assay Kit

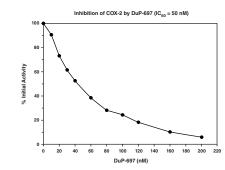
760151

- Measure cyclooxygenase (COX) activity in cell lysates, tissue homogenates, and purified enzyme preparations
- Measure the peroxidase component of the COX
- Assay 36 samples in duplicate
- Plate-based, colorimetric measurement (590 nm)
- Includes isozyme-specific inhibitors for distinguishing COX-2 activity from COX-1 activity



#### COX Colorimetric Inhibitor Screening Assay Kit 701050

- Screen for inhibitors of ovine cyclooxygenase (COX)-1 or human COX-2
- Measure the peroxidase component of the COX
- Assay 41 samples for COX-1 and 41 samples for COX-2 in duplicate
- Plate-based, colorimetric measurement (590 nm)



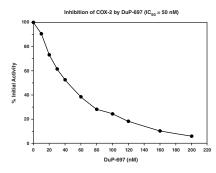
760111

COX (ovine) Inhibitor Screening Assay Kit

C-Reactive Protein (human) ELISA Kit

10011236

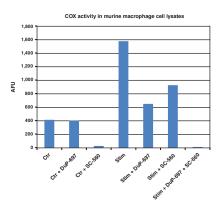
- Screen for inhibitors of ovine cyclooxygenase (COX)-1 or ovine COX-2
- Measure the peroxidase component of the COX
- Assay 41 samples for COX-1 and 41 samples for COX-2 in duplicate
- Plate-based, colorimetric measurement (590 nm)



#### **COX Fluorescent Activity Assay Kit**

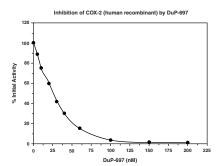
700200

- Measure cyclooxygenase (COX) activity in cell lysates, tissue homogenates, and purified enzyme preparations
- Measure the peroxidase component of the COX
- Assay 70 samples in duplicate
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)
- Includes isozyme-specific inhibitors for distinguishing COX-2 activity from COX-1 activity



#### COX Fluorescent Inhibitor Screening Assay Kit 700100

- Screen for inhibitors of ovine cyclooxygenase (COX)-1 or human COX-2
- Measure the peroxidase component of the COX
- Assay 46 samples for COX-1 and 46 samples for COX-2 in duplicate
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)

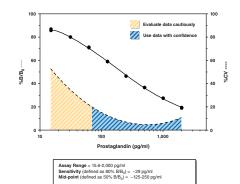


• Screen for inhibitors of ovine cyclooxygenase (COX)-1 or ovine COX-2

560101

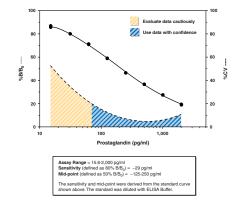
701070

- Directly measure PGF<sub>20</sub> produced in the COX reaction
- Assay 15 samples for COX-1 or COX-2 in duplicate • Plate-based, colorimetric measurement (405-420 nm)



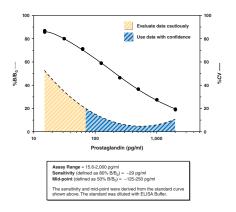
#### COX (ovine/human) Inhibitor Screening Assay Kit 560131

- Screen for inhibitors of ovine cyclooxygenase (COX)-1 or human COX-2
- Directly measure PGF produced in the COX reaction
- Assay 15 samples for COX-1 or COX-2 in duplicate
- Plate-based, colorimetric measurement (405-420 nm)



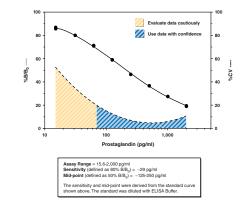
#### COX-1 (human) Inhibitor Screening Assay Kit

- Screen for inhibitors of human cyclooxygenase 1 (COX-1)
- Directly measure prostaglandin production
- Assay 16 samples in duplicate
- Plate-based, colorimetric measurement (405-420 nm)



#### COX-1 (ovine) Inhibitor Screening Assay Kit

- Screen for inhibitors of ovine cyclooxygenase 1 (COX-1)
- Directly measure prostaglandin production
- Assay 16 samples in duplicate
- Plate-based, colorimetric measurement (405-420 nm)

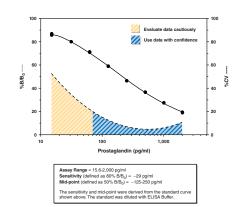


#### COX-2 (human) Inhibitor Screening Assay Kit

701080

701090

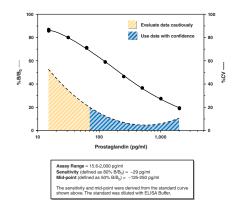
- Screen for inhibitors of human cyclooxygenase 2 (COX-2)
- Directly measure prostaglandin production
- Assay 16 samples in duplicate
- Plate-based, colorimetric measurement (405-420 nm)



#### COX-2 (ovine) Inhibitor Screening Assay Kit

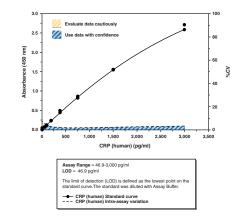
702000

- Screen for inhibitors of ovine cyclooxygenase 2 (COX-2)
- Directly measure prostaglandin production
- Assay 16 samples in duplicate
- Plate-based, colorimetric measurement (405-420 nm)



#### • Measure C-reactive protein (CRP) from human plasma without prior sample purification

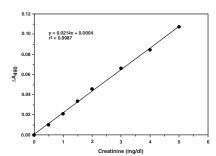
- Assay Range: 46.9-3,000 pg/ml with an LOD of 46.9 pg/ml
- Rapid assay; get results in under 3 hours



#### Creatinine (serum) Colorimetric Assay Kit

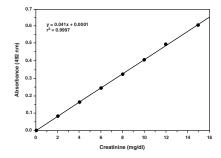
700460

- Measure creatinine levels in plasma and serum
- Assay 40 samples in duplicate
- Measure creatinine levels down to 0.1 mg/dl
- Plate-based, colorimetric measurement (490-500 nm)



#### Creatinine (urinary) Colorimetric Assay Kit

- Measure creatinine levels in urine
- Assay 40 samples in duplicate
- Measure creatinine levels down to 0.1 mg/dl
- Plate-based, colorimetric measurement (490-500 nm)



CREB (Phospho-Ser<sup>133</sup>) Transcription Factor **Assay Kit** 

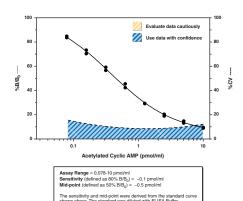
10009846

- A sensitive, non-radioactive method of detecting cAMP-responseelement-binding protein (CREB) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates

#### Cyclic AMP ELISA Kit

581001

- Measure acetylated or non-acetylated cyclic AMP (cAMP) in cell lysates, tissue homogenates, plasma or urine
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure cAMP levels down to 0.1 pmol/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm
- Contains acetic anhydride

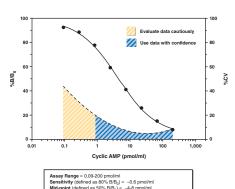


Also Available: Cyclic AMP ELISA Kit (without Acetic Anhydride) Item No. 581002

#### Cyclic AMP Select ELISA Kit

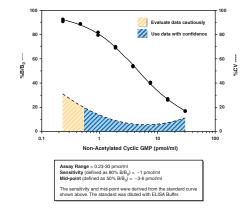
501040

- A sensitive assay for cyclic AMP (cAMP) in multiple sample matricies without the need for acetylation
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure cAMP levels down to 0.6 pmol/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



Cyclic GMP ELISA Kit

- Detect non-acetylated or acetylated cyclic GMP (cGMP) in cell lysates, tissue homogenates, plasma, or urine
- Measure non-acetylated cGMP levels down to 1 pmol/ml or acetylated cGMP levels down to 0.1 pmol/ml
- Assay 24 samples in triplicate or 36 samples in duplicate
- Contains acetic anhydride
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



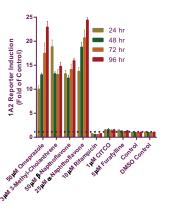
Also Available: Cyclic GMP ELISA Kit (without Acetic Anhydride) Item No. 581022

#### CYP1A1/2 Induction Reporter Assay Kit

600670

581021

- A 96-well plate coated with DNA encoding a CYP1A1/2 promoter-regulated reporter enzyme (SEAP)
- Cells grown on the complex will express the CYP1A1/2 promoter-regulated SEAP
- Binding of inducer-activated endogenous transcription factors to the CYP1A1/2 promoter will activate SEAP expression
- SEAP secreted into the supernatant cleaves a luminescence-based



CYP2B6 Induction Reporter Assay Kit

receptors, hCAR3 and HNF4α

substrate

• A 96-well plate coated with DNA encoding a CYP2B6

and the CYP2B6 promoter-regulated SEAP

promoter-regulated reporter enzyme (SEAP) and two nuclear

• CAR-mediated CYP2B6 inducers will activate SEAP expression

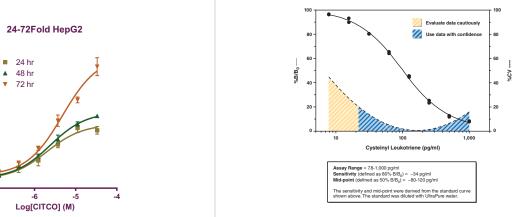
• Cells grown on the complex will express the two nuclear receptors

• SEAP secreted into the supernatant cleaves a luminescence-based

#### Cysteinyl Leukotriene ELISA Kit

500390

- Detect all cysteinyl leukotrienes (CysLTs) (LTC,, LTD,, and LTE,) in urine and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure CysLT levels down to 34 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

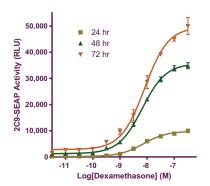


#### CYP2C9 Induction Reporter Assay Kit

601120

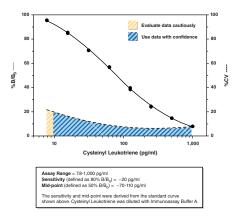
600680

- A 96-well plate coated with DNA encoding a CYP2C9 promoter-regulated reporter enzyme (SEAP) and two nuclear receptors, GRα and HNF4α
- Cells grown on the complex will express the two nuclear receptors and the CYP2C9 promoter-regulated SEAP
- GR-mediated CYP2C9 inducers will activate SEAP expression
- SEAP secreted into the supernatant cleaves a luminescence-based substrate



#### Cysteinyl Leukotriene Express ELISA Kit

- Detect all cysteinyl leukotrienes (CysLTs) (LTC,, LTD,, and LTE,) in urine, tissue culture medium, and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure CysLT levels down to 20 pg/ml
- Incubation: 2 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



• Assay 24 samples in triplicate or 36 samples in duplicate

Assay Range = 15.6-2,500 pg/ml Sensitivity (defined as 80% B/B<sub>o</sub>) = -70 pg/ml Mid-point (defined as 50% B/B<sub>o</sub>) = -200-300 pg/m

EP<sub>2</sub> Receptor (rat) Reporter Assay Kit

and the SEAP reporter

substrate

e sensitivity and mid-point were derived from the standa own above. The standard was diluted with ELISA Buffer.

• A 96-well plate coated with DNA encoding both the rat EP,

receptor and a cAMP response element-regulated SEAP

• Binding of EP, receptor agonists induces expression of SEAP

• Cells grown on the complex will express both the rat EP<sub>2</sub> receptor

SEAP secreted into the supernatant cleaves a luminescence-based

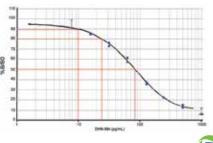
• Measure enterolactone levels down to 70 pg/ml

orbital shaker | Read: Colorimetric at 405-420 nm

• Measure enterolactone in plasma, urine, and other sample matrices

• Incubation: 18 hours at 4°C | Development: 90–120 minutes on an

- Measure 1,4-dihydroxynonane mercapturic acid (DHN-MA), a HNE
- metabolite, in urine samples • Assay 24 samples in triplicate or 36 samples in duplicate
- Incubation: Overnight
- Species Compatibility: Human | Mouse | Rat | Dog | Pig



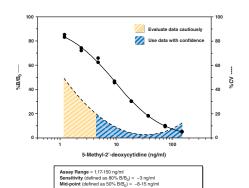


#### **DNA Methylation ELISA Kit**

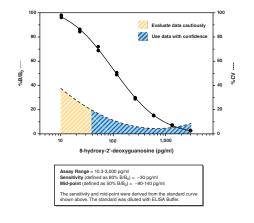
589324

501140

- Measure 5-methyl-2'-deoxy cytidine in urine, plasma, cell lysates, and tissue culture supernatants
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 5-methyl-2'-deoxy cytidine levels down to 3 ng/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



- Measure major oxidative damage markers, 8-hydroxy 2'-deoxyguanosine, 8-hydroxyguanosine, and 8-hydroxyguanine in urine, cell culture medium, cell lysates, tissue samples, saliva, and plasma/serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Assay Range: 10.3-3,000 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

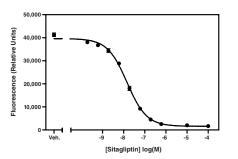


Also Available: 8-Hydroxyguanosine Monoclonal ELISA Kit Item No. 501130 (see page 42)

#### DPP (IV) Inhibitor Screening Assay Kit

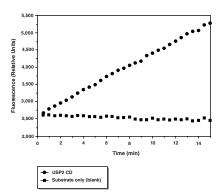
700210

- Screen for dipeptidyl peptidase IV (DPP (IV)) inhibitors
- Assay 46 samples in duplicate
- Includes positive control inhibitor
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### · Screen potential inhibitors and activators for activity against specific deubiquitinating enzymes (DUBs)

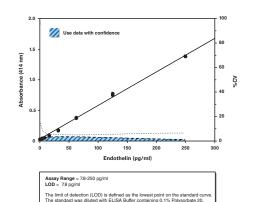
- · Assess performance of DUBs of interest with ubiquitin-AMC substrate allowing relevant kinetic data to be generated
- Evaluate performance of DUB activators and inhibitors enabling relevant kinetic data to be determined
- Optimize assays for specific DUBs to facilitate their use in high-throughput screening
- Demonstrate novel, putative DUBs have ubiquitin-AMC processing activity



#### **Endothelin ELISA Kit**

583151

- Measure endothelin in serum, plasma, urine, or cell culture media
- Assay 24 samples in triplicate or 36 samples in duplicate
- Assay Range: 7.8-250 pg/ml with an LOD of 7.8 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



#### EP, Receptor (rat) Activation Assay Kit (cAMP)

# 600410

600340

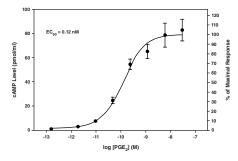
• Reverse transfection reporter assay to screen for EP, receptor agonists

log [PGE<sub>2</sub>] (M)

- Assay 18 samples in duplicate
- ELISA-based cAMP measurement

-- No AH6809 -- 60 μM AH6809

• Plate-based, colorimetric measurement (405-420 nm)



EP, Receptor (rat) Reporter Assay Kit

ERα (human) Reporter Assay Kit

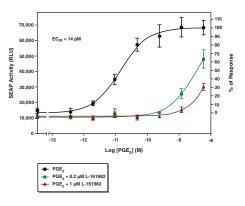
15737

ERRy (human) Reporter Assay Kit

FAAH Inhibitor Screening Assay Kit

10005196

- A 96-well plate coated with DNA encoding both the rat EP, receptor and a cAMP response element-regulated SEAP
- Cells grown on the complex will express both the rat EP, receptor and the SEAP reporter
- Binding of EP, receptor agonists induces expression of SEAP
- SEAP secreted into the supernatant cleaves a luminescence-based substrate

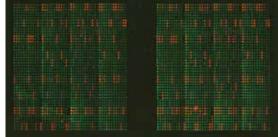


#### EpiTitan<sup>™</sup> Histone Peptide Array

17626

600350

- Array covers over 95 unique modifications on the four core histones
- Perform multiple experiments per array
- Determine the specificity of histone binding proteins, the selectivity of histone antibodies and identify substrates for histone-modifying



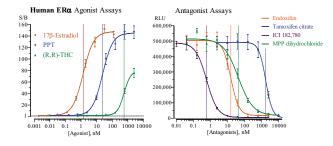
#### ER (human) Reporter Assays Panel

15739

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

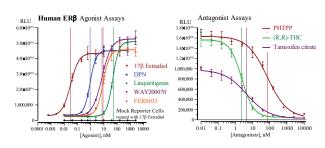




#### ERβ (human) Reporter Assay Kit

15738

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

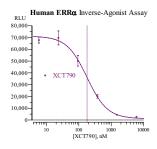




#### ERRα (human) Reporter Assay Kit

15760

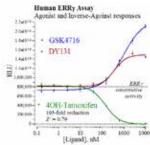
- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details





#### • A nuclear receptor cell-based reporter assay

• See www.caymanchem.com for more details



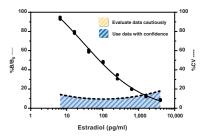


#### **Estradiol ELISA Kit**

582251

16524

- Measure estradiol (17β-estradiol) in cell culture medium, plasma, and serum
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure estradiol levels down to 15 pg/ml
- Incubation: 1 hour | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm

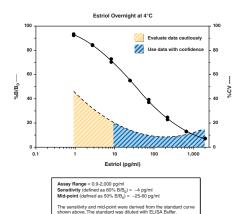


ssay Range = 6.6-4,000 pg/ml ensitivity (defined as 80% B/B<sub>0</sub>) = ~15 pg/ml

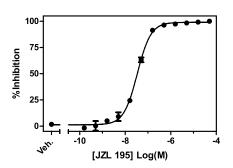
#### **Estriol ELISA Kit**

582281

- Measure estriol, a major estrogen produced during pregnancy, in plasma and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure estriol levels down to 4 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



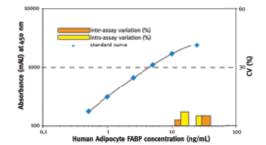
- Screen for inhibitors of fatty acid amide hydrolase (FAAH)
- Assay 45 samples in duplicate
- Includes JZL 195 positive control
- Plate-based, fluorometric measurement (ex 340-360 nm, em 450-465 nm)



#### FABP4 (human) EIA Kit

10007614

- Measure fatty acid binding protein 4 (FABP4) in human serum, plasma, tissue extract, and culture supernatant
- Assay 41 samples in duplicate
- Incubations: 2 x 1 hour | 30 minutes



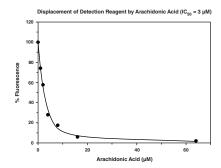


10010231

#### FABP4 Inhibitor/Ligand Screening Assay Kit

• Screen for fatty acid binding protein 4 (FABP4) ligands

- Includes arachidonic acid positive control
- Plate-based, fluorometric measurement (ex 370 nm, em 475 nm)



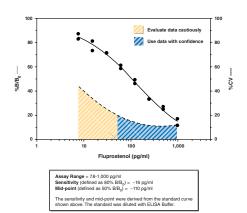
Fluprostenol ELISA Kit

#### FXR (human) Reporter Assay Kit

#### 10006306

#### 10006307

- Measure the free acid form of fluprostenol
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure fluprostenol levels down to 16 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

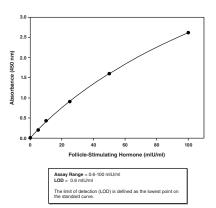


#### Follicle-Stimulating Hormone ELISA Kit

500710

516761

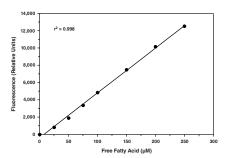
- Measure follicle-stimulating hormone (FSH) in human serum
- Assay 27 samples in triplicate or 41 samples in duplicate
- Assay Range: 0.6-100 mIU/mI with an LOD of 0.6 mIU/mI
- Rapid assay; get results in under 2 hours



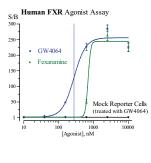
#### Free Fatty Acid Fluorometric Assay Kit

700310

- Measure free fatty acids (FFAs) in plasma, serum, and urine
- Assav 20 samples in duplicate
- Measure free fatty acids from 25 to 250  $\mu M$
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



# **MINDIGO**

#### GFAP (human) EIA Kit

10007621

15741

- Measure glial fibrillary acidic protein (GFAP) in human serum, plasma, CSF, and tissue culture supernatant samples
- Assay 40 samples in duplicate
- Incubations: 2 hours | 2 x 1 hour



#### Acylated Ghrelin (dog) Easy Sampling EIA Kit

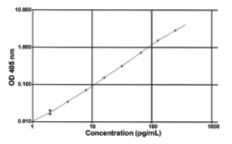
501210

- Measure acetylated ghrelin in dog blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubations: 1 hour | Overnight
- Samples collected on any protease inhibitor are suitable with this kit, including frozen samples



#### Acylated Ghrelin (human) Easy Sampling EIA Kit

- Measure acetylated ghrelin in human blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubation: 2 x 2 hours
- Samples collected on any protease inhibitor are suitable with this kit, including frozen samples



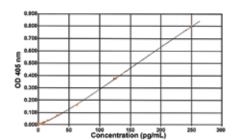


# • Measure acetylated ghrelin in human blood, plasma, or serum

• Assay 24 samples in triplicate or 36 samples in duplicate

Acylated Ghrelin (human) Express EIA Kit

- **Incubation:** 3 hours
- Sampling tubes with PHMB available to store, preserve, and protect samples from protein degradation (see Item No. 10949)

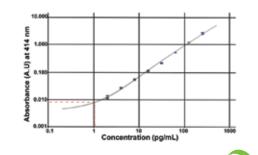


#### Acylated Ghrelin (mouse, rat) Easy Sampling EIA Kit

501170

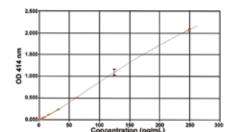
SPI bio beptin

- Measure acetylated ghrelin in mouse or rat blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- **Incubation:** 2 x 2 hours
- Samples collected on any protease inhibitor are suitable with this kit, including frozen samples



# Acylated Ghrelin (mouse, rat) Express EIA Kit

- Measure acetylated ghrelin in mouse or rat blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubation: 3 hours
- Sampling tubes with PHMB available to store, preserve, and protect samples from protein degradation (see Item No. 10949)





#### Acylated Ghrelin (pig) EIA Kit

501220

- Measure acetylated ghrelin in pig blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubation: Overnight
- Sampling tubes with PHMB available to store, preserve, and protect samples from protein degradation (see Item No. 10949)





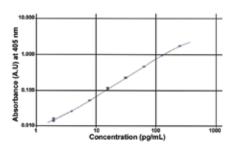
#### Unacylated Ghrelin (dog) Easy Sampling EIA Kit

- Measure unacetylated ghrelin in dog blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubations: 1 hour | Overnight
- Samples collected on any protease inhibitor are suitable with this kit, including frozen samples





- 501190
- Measure unacetylated ghrelin in human blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- **Incubation:** 2 x 2 hours
- Samples collected on any protease inhibitor are suitable with this kit, including frozen samples

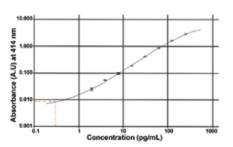




#### Unacylated Ghrelin (mouse, rat) Easy Sampling EIA Kit

501180

- Measure unacetylated ghrelin in mouse or rat blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- **Incubation:** 2 x 2 hours
- Samples collected on any protease inhibitor are suitable with this kit, including frozen samples





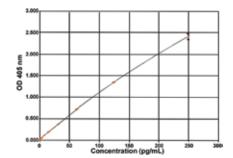
#### Unacylated Ghrelin (pig) EIA Kit

501230

- Measure unacetylated ghrelin in pig blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubation: Overnight
- Sampling tubes with PHMB available to store, preserve, and protect samples from protein degradation (see Item No. 10949)



- Measure unacetylated ghrelin in human blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubation: 3 hours
- Collection tubes with PHMB available for easy sample collection (see Item No. 10949)

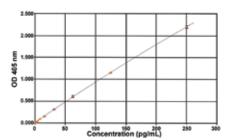




#### Unacylated Ghrelin (mouse, rat) **Express EIA Kit**

10008953

- Measure unacetylated ghrelin in mouse or rat blood, plasma, or serum samples
- Assay 24 samples in triplicate or 36 samples in duplicate
- Incubation: 3 hours
- Sampling tubes with PHMB available to store, preserve, and protect samples from protein degradation (see Item No. 10949)

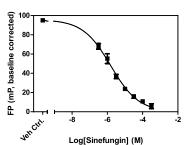




#### GLP SAM-Screener<sup>™</sup> Assay Kit

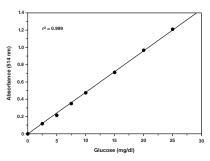
600570

- Homogeneous, fluorescence polarization assay to screen for SAMbinding site inhibitors
- Designed for high-throughput screening
- Plate-based, fluorescence polarization measurement (ex 575 nm,



#### • Measure glucose in plasma, serum, and urine

- Assay 40 samples in duplicate
- Assay Range: 2.5-25 mg/dl
- Plate-based, colorimetric measurement (500-520 nm)



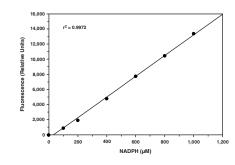
#### Glucose-6-Phosphate Dehydrogenase **Activity Assay Kit**

700300

700750

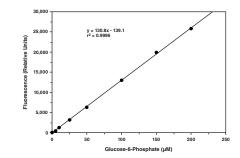
10009582

- Measure glucose-6-phosphate dehydrogenase (G6PDH) activity in cell lysates and tissue homogenates
- Assay 20 samples in duplicate
- Measure G6PDH activity down to 5 U/ml
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)

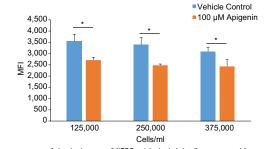


#### Glucose-6-Phosphate Fluorometric Assay Kit

- Measure glucose-6-phosphate (G6P) in cell lysates and tissue homogenates
- Assay 20 samples in duplicate
- Assay Range: 5-200 μM
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



- Measure glucose uptake by cultured cells
- Employ 2-NBDG, a fluorescent deoxyglucose analog • An inhibitor of glucose transport is included as a control
- Convenient tool for studying modulators of glucose uptake



Apigenin decreases 2-NBDG uptake in Jurkat cells as measured by flow cytometry.

#### Glutaredoxin Fluorescent Activity Assay Kit

• Convenient 96-well plate format

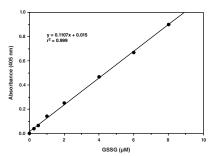
- Measure glutathione reductase activity in cell lysates and tissue homogenates
- Active human glutaredoxin 1 provided with kit



11536

#### Glutathione Assay Kit

- Measure total, oxidized glutathione (GSSG), and/or reduced glutathione (GSH) in cell lysates, tissue homogenates, plasma and erythrocyte lysates, and serum
- Assay 40 samples in duplicate
- Assay Range: 0.25-8  $\mu$ M (GSSG) or 0.5-16  $\mu$ M (GSH)



Glutathione Cell-Based Detection Kit (Blue Fluorescence)

(GSH) probe

cell lysates

600360

- Measure glutathione S-transferase (GST) activity in plasma, tissue homogenates, and cell lysates • Utilize monochlorobimane (MCB), a highly fluorescent glutathione
  - Assay 22 samples in duplicate
  - Measure GST activity down to 24 U/ml

S-Glutathionylated Protein Detection Kit

• Reagents provided to test up to 30 samples

Glycerol Cell-Based Assay Kit

produces a brilliant purple product

acid cycling rate

extracellular glycerol

• Direct visualization of S-glutathionylated proteins in whole

(permeabilized) cells by flow cytometry and microscopy • Visualization using colorimetric or fluorescence detection

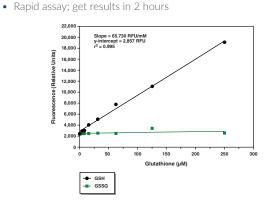
Mouse monocytes – Fluorescence microscopy. Cells were stained with avidin-FITC using protoco supplied. Panel A: No reduction reagent. Panel B: Full protocol reveals S-glutathionylated proteins. Panel C No free-thiol Blocking Reagent reveals labeling of all accessible protein thiols.

• Detect extracellular glycerol as a measure of the triglyceride/fatty

• Measure glycerol by a coupled enzymatic reaction system that

• Includes a standard curve for accurate quantification of

• Plate-based, colorimetric measurement (340 nm)

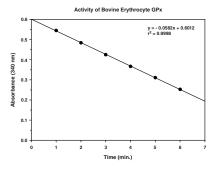


• Includes a standard curve for accurate quantification of GSH from

#### Glutathione Peroxidase Assay Kit

703102

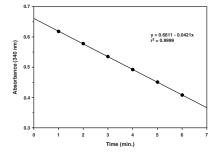
- Measure glutathione peroxidase (GPx) in plasma, erythrocyte lysate, tissue homogenates, and cell lysates
- Assay 30 samples in triplicate or 45 samples in duplicate
- Measure GPx activity down to 50 U/ml



#### Glutathione Reductase Assay Kit

703202

- Measure glutathione reductase (GR) activity in plasma, cell lysates, and tissue homogenates
- · Assay 45 samples in duplicate
- Measure GR activity down to 20 U/ml
- Plate-based colorimetric measurement (340 nm)



#### Glutathione S-Transferase Assay Kit

703302

10010721

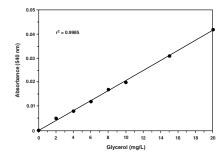
10011725

• Measure glycerol in plasma and serum

- Assay 40 samples in duplicate
- Assay Range: 2-20 mg/L

Glycerol Colorimetric Assay Kit

• Plate-based, colorimetric measurement (530-550 nm)

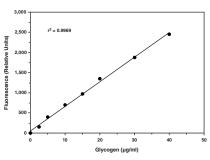


#### Glycogen Assay Kit

700480

10010755

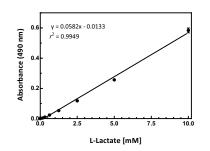
- Measure glycogen from tissue samples
- · Assay 20 samples in duplicate
- Assay Range: 2.5-40 μg/ml
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



# Glycolysis Cell-Based Assay Kit

600450

- Detect L-lactate, the end product of glycolysis
- Can be adapted to high-throughput screening
- Measure L-lactate in cell culture supernatant down to 156 µM



#### Glyco-SPOT DNA Repair Assay Kit

• Profile glycosylase/AP endonuclease activities

- Media Application: Pure Enzyme | Cell Nuclear Extracts | Whole Cell Extracts | Tissue Extracts
- HTS friendly/fluorescent micro array
- Species Compatibility: Human | Mouse | Rat | Fish | Bacteria



501280

#### Goat Anti-Human Fluorophore Selection Kit

16541

- Comprehensive: Columbia Biosciences' fluorophore selection kits include the secondary antibody individually conjugated to seven unique fluorophores
- Efficient: The ability to evaluate the various labeled conjugates side-by-side allows scientists to rapidly optimize their reagents
- Cost-effective: Order one kit and get seven conjugated secondary antibodies. The fluorophore selection kit reduces costs associated with purchasing and shipping full-size antibody conjugates.





#### Goat Anti-Mouse Fluorophore Selection Kit

- Comprehensive: Columbia Biosciences' fluorophore selection kits include the secondary antibody individually conjugated to seven unique fluorophores
- Efficient: The ability to evaluate the various labeled conjugates side-by-side allows scientists to rapidly optimize their reagents
- Cost-effective: Order one kit and get seven conjugated secondary antibodies. The fluorophore selection kit reduces costs associated with purchasing and shipping full-size antibody conjugates.





- Comprehensive: Columbia Biosciences' fluorophore selection kits include the secondary antibody individually conjugated to seven unique fluorophores
- Efficient: The ability to evaluate the various labeled conjugates side-by-side allows scientists to rapidly optimize their reagents
- Cost-effective: Order one kit and get seven conjugated secondary antibodies. The fluorophore selection kit reduces costs associated with purchasing and shipping full-size antibody conjugates.



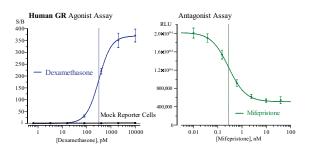


#### GR (human) Reporter Assay Kit

15733

16543

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

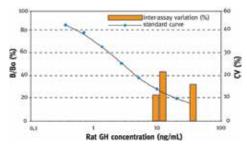




#### Growth Hormone (rat) EIA Kit

589601

- Measure growth hormone (GH) in rat plasma, and cell culture supernatant samples
- Assay 33 samples in duplicate
- Incubation: 2 x Overnight
- Calibrated against the reference preparation NIDDK standard rGH-RP2



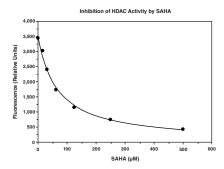


- Screen for inhibitors of pCAF
- Assay 46 samples in duplicate
- Plate-based, fluorometric measurement (ex 360-390 nm, em 450-470 nm)

#### **HDAC Cell-Based Activity Assay Kit**

600150

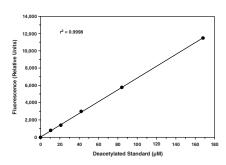
- Screen for modulators of histone deacetylase (HDAC) activity in whole cells
- Includes trichostatin A as a positive control
- Plate-based, fluorometric measurement (ex 340-360 nm, em 440-460 nm)



#### **HDAC Fluorometric Activity Assay Kit**

10011563

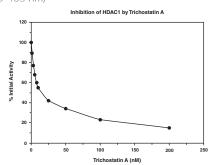
- Measure class I and class II histone deacetylase (HDAC) activity in nuclear extracts
- Assay 20 samples in duplicate
- Measure HDAC activity down to 7 U/ml
- Plate-based, fluorometric measurement (ex 340-360 nm, em 440-465 nm)



#### **HDAC1** Inhibitor Screening Assay Kit

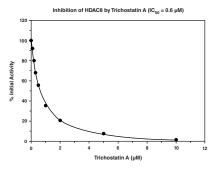
10011564

- Screen for inhibitors of histone deacetylase 1 (HDAC1)
- Assay 45 samples in duplicate
- Includes trichostatin A as a positive control
- Plate-based, fluorometric measurement (ex 340-360 nm, em 440-465 nm)



#### **HDAC8** Inhibitor Screening Assay Kit

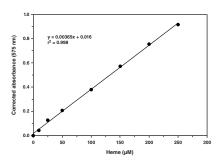
- Screen for inhibitors of histone deacetylase 8 (HDAC8)
- Assay 45 samples in duplicate
- Includes trichostatin A as a positive control
- Plate-based, fluorometric measurement (ex 340-360 nm, em 440-465 nm)



#### Hemoglobin Colorimetric Assay Kit

700540

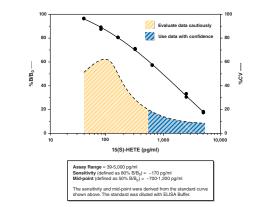
- Measure total hemoglobin concentration in blood, tissue homogenates, and cell lysates
- Assay 40 samples in duplicate
- Assay Range: 0.016-0.4 g/dl
- Plate-based, colorimetric measurement (560-590 nm)



#### 15(S)-HETE ELISA Kit

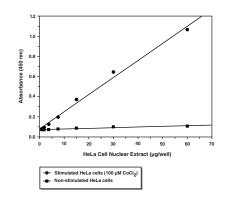
534721

- Detect 15(S)-HETE, a product of 15-lipoxygenase treatment of arachidonic acid
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 15(S)-HETE levels down to 170 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



### HIF-1α Transcription Factor Assay Kit

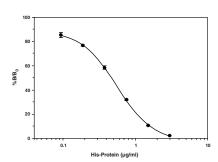
- A sensitive, non-radioactive method for detecting hypoxia inducible factor- $1\alpha$  (HIF- $1\alpha$ ) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays (EMSAs)
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



#### His-Tag Detection ELISA Kit

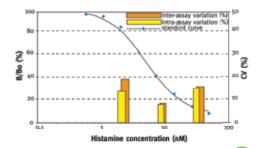
10012445

- A semi-quantitative screening assay to detect hexahistidine-tagged (His-tagged) proteins from cell lysates or affinity column fractions
- An alternative to SDS-PAGE for testing affinity column fractions • Assay 26 samples in triplicate or 39 samples in duplicate
- Rapid assay; get results in under 3 hours



#### Histamine EIA Kit

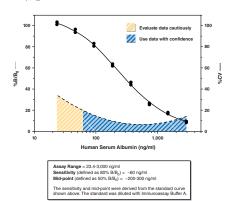
- Measure histamine in plasma, serum, bronchoalveolar lavage, tissue extracts, and cell culture supernatant samples
- Assay 33 samples in duplicate
- Incubation: Overnight







- Measure the presence of albumin in human urine, an indicator of renal function
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure albumin levels down to 60 ng/ml
- Rapid assay; get results in under 4 hours

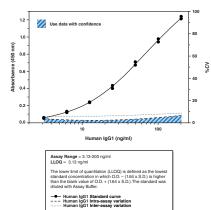


#### Human Therapeutic IgG1 ELISA Kit

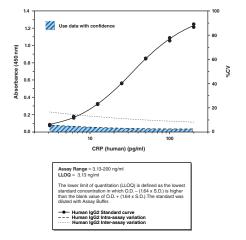
500910

500340

- Specifically detect human IgG1 antibodies in monkey plasma or serum
- Not cross reactive to rhesus or cynomolgus immunoglobulins
- Useful for pharmacokinetic analysis of therapeutic human IgG1 antibodies in monkeys
- Useful for detecting human IgG1 in mouse or rat plasma or serum



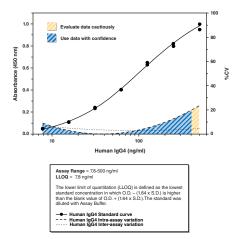
- Specifically detect human IgG2 antibodies in monkey plasma or serum
- Not cross reactive to rhesus or cynomolgus immunoglobulins
- Useful for pharmacokinetic analysis of therapeutic human IgG2 antibodies in monkeys
- Useful for detecting human IgG2 in mouse or rat plasma or serum



#### Human Therapeutic IgG4 ELISA Kit

500920

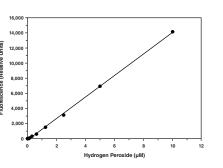
- Specifically detect human IgG4 antibodies in monkey plasma or serum
- Not cross reactive to rhesus or cynomolgus immunoglobulins
- Useful for pharmacokinetic analysis of therapeutic human IgG4 antibodies in monkeys
- Useful for detecting human IgG4 in mouse or rat plasma or serum



#### Hydrogen Peroxide Cell-Based Assay Kit

600050

- A simple fluorometric assay for the quantification of extracellular H<sub>2</sub>O<sub>2</sub>
- Utilize ADHP, a sensitive and stable probe for H<sub>2</sub>O<sub>2</sub>
- Catalase, an H<sub>2</sub>O<sub>2</sub> scavenger, is included as a control for specificity
- Rapid assay; results in under 1 hour

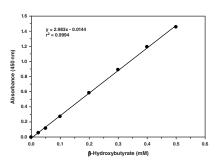


# Colorimetric Assay Kit

- Measure  $\beta$ -hydroxybutyrate ( $\beta$ -HB) from plasma, serum, urine, cell lysates, and tissue homogenates
- Assay 40 samples in duplicate

β-Hydroxybutyrate (Ketone Body)

- Assay Range: 25-500 μM
- Plate-based colorimetric measurement (445-455 nm)

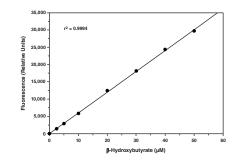


#### β-Hydroxybutyrate (Ketone Body) Fluorometric Assay Kit

700740

700190

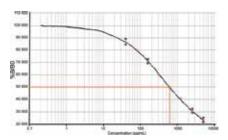
- Measure β-hydroxybutyrate (β-HB) from plasma, serum, urine, cell lysates, and tissue homogenates
- Assay 20 samples in duplicate
- Assay Range: 2.5-50 μM
- Plate-based fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### 20-Hydroxyecdysone EIA Kit

501390

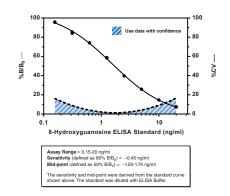
• Exclusive assay to measure 20-hydroxyecdysone in insects, mollusks, other arthropods, extracts, other biological samples





# 8-Hydroxyguanosine Monoclonal ELISA Kit

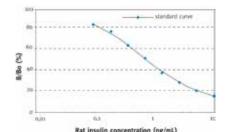
- Measure major oxidative damage markers, 8-hydroxy 2'-deoxyguanosine and 8-hydroxyguanosine in urine and other sample matrices as verified by MS analysis
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 8-hydroxyguanosine levels down to 0.45 ng/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



Also Available: DNA/RNA Oxidative Damage ELISA Kit Item No. 589320 (see page 29)

#### Insulin (rat) EIA Kit

- Measure insulin in human, mouse, or rat plasma, serum, and cell culture supernatant samples
- Assay 33 samples in duplicate
- Incubation: Overnight
- Insulinase Inhibitor Cocktail (Item No. 18429) available for easy and stable sample collection





Interleukin-1α (human) ELISA Kit

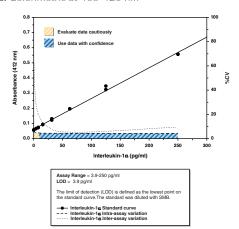
#### Interleukin-6 (human) ELISA Kit

#### 501030 8-Isoprostane ELISA Kit

#### JMJD2A Tudor Domains TR-FRET Assay Kit

700960

- Measure interleukin-1α (IL-1α) from human plasma, serum, synovial fluid, and other sample matrices
- Assay Range: 3.9-250 pg/ml with an LOD of 3.9 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm

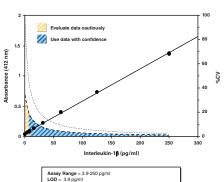


#### Interleukin-1ß (human) ELISA Kit

583311

583301

- Measure interleukin-1β (IL-1β) from human plasma, serum, synovial fluid, and other sample matrices
- Assay Range: 3.9-250 pg/ml with an LOD of 3.9 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm

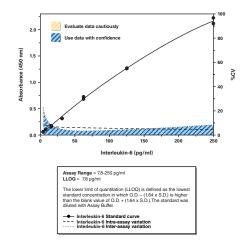




• Measure interleukin-6 (IL-6) from human plasma, serum, synovial fluid, and other sample matrices

• Assay Range: 7.8-250 pg/ml with an LLOQ of 7.8 pg/ml

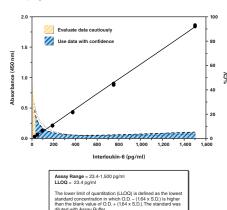
• Rapid assay; get results in under 3 hours



#### Interleukin-6 (mouse) ELISA Kit

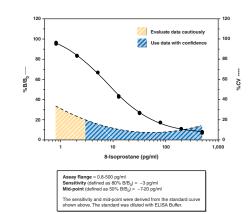
583371

- Measure interleukin-6 (IL-6) from mouse plasma, serum, and other sample matrices
- Assay Range: 23.4-1,500 pg/ml with an LLOQ of 23.4 pg/ml
- Rapid assay; get results in under 3 hours



#### • Measure 8-isoprostane, a biomarker of oxidative stress and antioxidant deficiency

- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 8-isoprostane levels down to 3 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

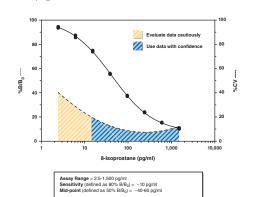


#### 8-Isoprostane Express ELISA Kit

516360

516351

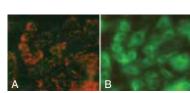
- Measure 8-isoprostane, a biomarker of oxidative stress and antioxidant deficiency
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 8-isoprostane levels down to 10 pg/ml
- Rapid assay; get results in under 4 hours



#### JC-1 Mitochondrial Membrane Potential **Assay Kit**

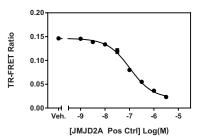
10009172

- Measure mitochondrial membrane potential as an indicator of cell health
- Utilize JC-1, a fluorescent, lipophilic, cationic dye
- Red fluorescence indicates healthy mitochondria and green fluorescence indicates mitochondria in poor health
- Staining in less than 1 hour



H9C2 cells stained with JC-1 followed by FCCP treatment Panel A: Cells prior to FCCP treatment shows JC-1 J-aggregates (red) accumulated in the mitochondria. Panel B: Cells subsequent to FCCP treatment shows diffuse JC-1 J-monomers (green) in the mitochondria.

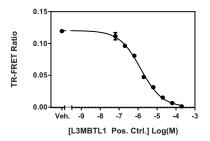
- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### L3MBTL1 MBT Domains TR-FRET Assay Kit

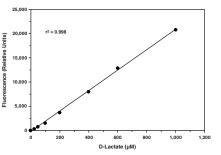
601030

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### **D-Lactate Assay Kit**

- Measure D-lactate in biological samples such as serum, plasma, whole blood, urine, saliva, and cell culture
- Assav 40 samples in duplicate
- Assay Range: 25 μM-1 mM
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



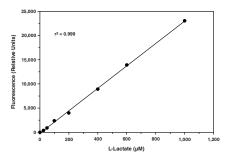
L-Lactate Assay Kit

LDH Cytotoxicity Assay Kit

601170

10007609 Leukotriene B, Express ELISA Kit

- Measure L-lactate in biological samples such as serum, plasma, whole blood, urine, saliva, and cell culture
- Assay 40 samples in duplicate
- Assay Range: 25 μM-1 mM
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)

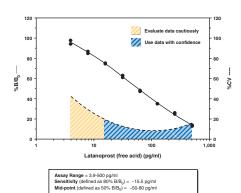


#### Latanoprost ELISA Kit

516811

700510

- Measure the free acid form of latanoprost
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure latanoprost levels down to 15.5 pg/ml
- Incubation: 2 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



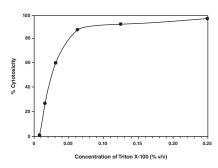
#### LC3 Interact Kit

15977

- Capture and detect LC3/GABARAP binding and associated proteins from biological samples
- Determine LC3/GABARAP selectivity of specific interacting proteins in vitro
- Identify LC3/GABARAP binding and associated proteins by western blotting or proteomic analysis
- Investigate the role of specific LC3 isoforms in different autophagy processes

• A colorimetric, non-radioactive assay for assessing cell damage

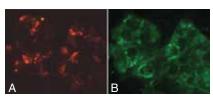
- Lactate dehydrogenase (LDH) release from the cytosol into the medium catalyzes a color-forming reaction
- Detect lysis of 10,000-200,000 cells per well
- Alternative to 51Cr-release assays



#### LDL Uptake Cell-Based Assay Kit

10011125

- Measure low-density lipoprotein (LDL) uptake in cultured cells
- Designed for fluorescence microscopy
- Measure LDL at ex 540 nm. em 570 nm
- Measure LDL-receptor at ex 485 nm, em 535 nm

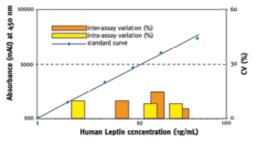


EGCG treatment of HepG2 cells stimulates uptake of LDL-DyLight™ 550. Panel A: LDL-DyLight™ 550 uptake is shown in red. Panel B: LDL

#### Leptin (human) EIA Kit

500010

- Measure leptin in human serum, plasma, and culture supernatant samples
- Assay 41 samples in duplicate
- Incubation: 2 x 1 hour

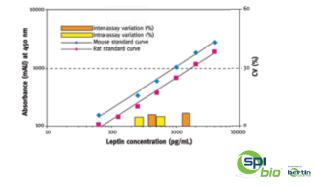




#### • Measure leptin in mouse or rat serum, plasma samples, and tissue culture supernatant

- Assay 40 samples in duplicate
- Incubation: 2 x 1 hour

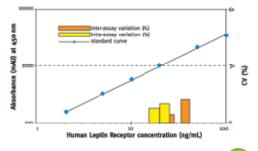
Leptin (mouse/rat) EIA Kit



#### Leptin Receptor (human) EIA Kit

10007608

- Measure leptin in human serum, plasma and culture supernatant
- Assay 41 samples in duplicate
- Incubation: 2 x 1 hour

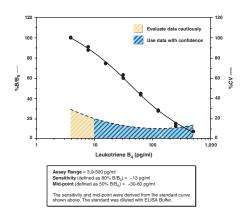


# SPI BOTTON

#### Leukotriene B<sub>4</sub> ELISA Kit

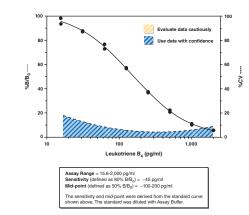
520111

- Measure leukotriene B<sub>4</sub> (LTB<sub>4</sub>) in plasma and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure LTB, levels down to 13 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



#### • Measure leukotriene B, (LTB,) in plasma and other sample matrices

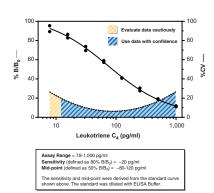
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure LTB, levels down to 45 pg/ml
- Rapid assay; get results in under 5 hours



#### Leukotriene C₁ ELISA Kit

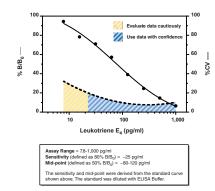
501070

- Measure leukotriene C, (LTC,) in tissue culture medium
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure LTC<sub>4</sub> levels down to 20 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



#### Leukotriene E₄ ELISA Kit

- Measure leukotriene  $E_4$  (LTE $_4$ ) in urine, with no purification required
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure LTE, levels down to 25 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



Lipase Activity Assay Kit

#### Lipid Hydroperoxide (LPO) Assay Kit (96 well)

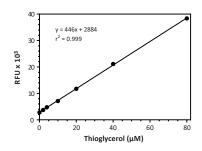
705003

Luteinizing Hormone ELISA Kit

Malachite Green Phosphate Assay Kit

600180

- Measure lipase activity in plasma, serum, tissue homogenates, and cell lysates
- Assay 20 samples in duplicate
- Measure lipase activity down to 0.1 U/ml
- Plate-based, fluorometric measurement (ex 380-390 nm, em 510-520 nm)

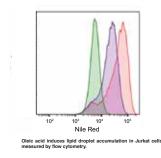


#### Lipid Droplets Fluorescence Assay Kit

#### 500001

700640

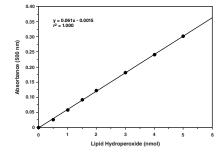
- Detect lipid droplets, cellular organelles that are also referred to as lipid bodies, oil bodies, or adiposomes
- Utilize Nile Red, a green fluorescent probe specific for lipid droplets
- Assess by fluorescence microscopy, flow cytometry, or plate reader
- Includes oleic acid as a positive control



#### Lipid Hydroperoxide (LPO) Assay Kit

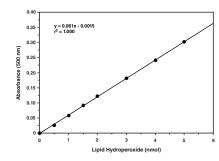
705002

- · Measure LPOs in samples
- Assay 40 samples in duplicate
- Assay Range: 0.25-5 nmol
- Plate-based, colorimetric measurement (500 nm)
- Requires reusable glass plate



• Measure LPOs in samples

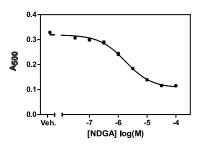
- Assay 40 samples in duplicate
- Assay Range: 0.25-5 nmol
- Plate-based, colorimetric measurement (500 nm)
- Includes reusable glass plate



#### Lipoxygenase Inhibitor Screening Assay Kit

#### 760700

- Screen for inhibitors of purified 5-, 12-, or 15-lipoxygenase (LO)
- Includes 15-LO
- Measure the hydroperoxides generated from LOs
- Plate-based, colorimetric measurement (490-500 nm)



#### Live Cell Multiplex Assay Kit

#### 15766

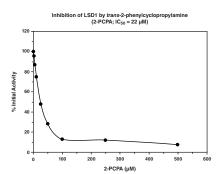
- · A fluorescence-based cell viability assay
- See www.caymanchem.com for more details



#### LSD1 Inhibitor Screening Assay Kit

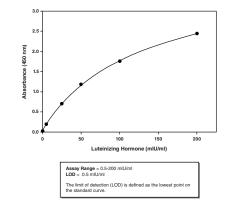
#### 700120

- Screen for inhibitors of lysine-specific demethylase 1 (LSD1)
- Measure the H<sub>2</sub>O<sub>2</sub> produced during demethylation
- Assay 45 samples in duplicate
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### • Measure luteinizing hormone (LH) in human serum

- Assay 27 samples in triplicate or 41 samples in duplicate • Assay Range: 0.5-200 mIU/mI with an LOD of 0.5 mIU/mI
- Rapid assay; get results in under 2 hours



#### LXR (human) Reporter Assays Panel

#### 15736

500720

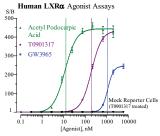
- · A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



#### LXRα (human) Reporter Assay Kit

#### 15735

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

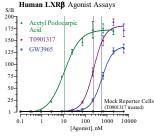


# **MINDIGO**

#### LXRβ (human) Reporter Assay Kit

#### 15734

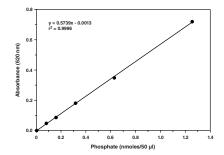
- A nuclear receptor cell-based reporter assay
- · See www.caymanchem.com for more details





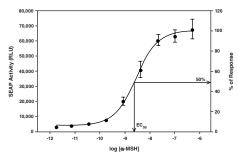
#### • Measure inorganic free phosphate in aqueous solutions

- Assay 42 samples in duplicate
- Assay Range: 1.6-25 μM
- Plate-based, colorimetric measurement (620-640 nm)



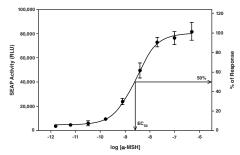
#### Melanocortin-3 Receptor Reporter Assay Kit

- Reverse transfection reporter assay to screen for melanocortin-3 receptor (MC3R) agonists
- Assay 46 samples in duplicate
- Includes α-MSH positive control
- Plate-based, luminometric measurement

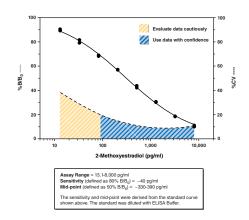


#### Melanocortin-4 Receptor Reporter Assay Kit 600190

- Reverse transfection reporter assay to screen for melanocortin-4 receptor (MC4R) agonists
- Assay 46 samples in duplicate
- Includes α-MSH positive control
- Plate-based, luminometric measurement



- Detect 2-methoxyestradiol (2-ME2)
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 2-ME2 levels down to 40 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

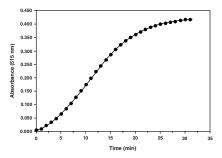


#### Methyltransferase Colorimetric Assay Kit

700140

582261

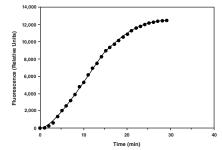
- Continuous assay to monitor SAM-dependent methyltransferases
- Assay 45 samples in duplicate
- Plate-based, colorimetric measurement (500-520 nm)



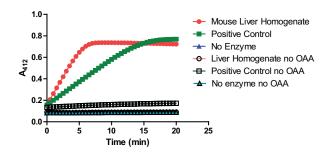
#### Methyltransferase Fluorometric Assay Kit

700150

- Continuous assay to monitor SAM-dependent methyltransferases
- Assay 36 samples in duplicate
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



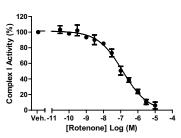
- Measure citrate synthase activity in isolated mitochondria, tissue and cell homogenate
- Assess mitochondrial content
- Suitable for high-throughput screening
- Plate-based, colorimetric measurement (412 nm)
- MitoCheck® Mitochondrial (Tissue) Isolation Kit (Item No. 701010) available to aid in mitochondrial tissue isolation



#### MitoCheck® Complex I Activity Assay Kit

700930

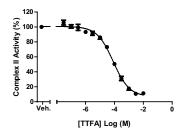
- Measure complex I activity in isolated mitochondria
- Useful for screening compounds for complex I inhibition
- Suitable for high-throughput screening
- Plate-based, colorimetric measurement (340 nm)
- MitoCheck® Mitochondrial (Tissue) Isolation Kit (Item No. 701010) available to aid in mitochondrial tissue isolation



#### MitoCheck® Complex II Activity Assay Kit

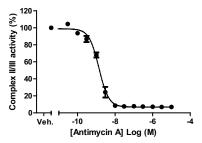
700940

- Measure complex II activity in isolated mitochondria
- Useful for screening compounds for complex II inhibition
- Suitable for high-throughput screening
- Plate-based, colorimetric measurement (600 nm)
- MitoCheck® Mitochondrial (Tissue) Isolation Kit (Item No. 701010) available to aid in mitochondrial tissue isolation



#### • Measure complex II/III activity in isolated mitochondria

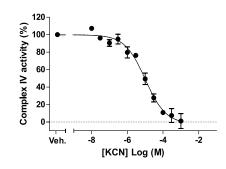
- Useful for screening compounds for complex II/III inhibition
- Suitable for high-throughput screening
- Plate-based, colorimetric measurement (550 nm)
- MitoCheck® Mitochondrial (Tissue) Isolation Kit (Item No. 701010) available to aid in mitochondrial tissue isolation



#### MitoCheck® Complex IV Activity Assay Kit

700990

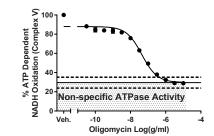
- Measure complex IV activity in isolated mitochondria
- Useful for screening compounds for complex IV inhibition
- Suitable for high-throughput screening
- Plate-based, colorimetric measurement (550 nm)
- MitoCheck® Mitochondrial (Tissue) Isolation Kit (Item No. 701010) available to aid in mitochondrial tissue isolation



#### MitoCheck® Complex V Activity Assay Kit

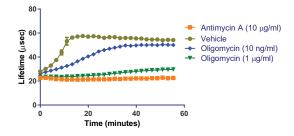
701000

- Measure complex V activity in isolated mitochondria
- Useful for screening compounds for complex V inhibition
- Suitable for high-throughput screening
- Plate-based, colorimetric measurement (340 nm)
- MitoCheck® Mitochondrial (Tissue) Isolation Kit (Item No. 701010) available to aid in mitochondrial tissue isolation



#### MitoCheck® Mitochondrial OCR Assay Kit

- Measure oxygen consumption rate (OCR) directly in freshly isolated mitochondria
- Useful for screening mitochondrial inhibitors and uncouplers
- Suitable for high-throughput screening
- Plate-based, fluorometric or time-resolved, fluorometric measurement (ex 380 nm, em 650 nm)
- MitoCheck® Mitochondrial (Tissue) Isolation Kit (Item No. 701010) available to aid in mitochondrial tissue isolation

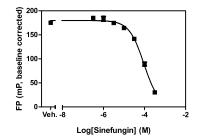


Sample Data. Oxygen consumption is measured in isolated mouse iver mitochondria (0.12 mg/ml) in the presence the complex V hnibitor oligonymic in 1 ug/ml, 10 ng/ml), the complex III inhibitor antimycin A (10 µM) and vehicle control. The difference in oxygen consumption rates in the absence and presence of oligomycin indicate that these mitochondria are coupled.

#### MLL1 SAM-Screener™ Assay Kit

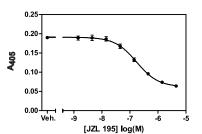
600580

- Homogeneous, fluorescence polarization assay to screen for SAMbinding site inhibitors
- Designed for high-throughput screening
- Plate-based, fluorescence polarization measurement (ex 575 nm, em 620 nm)



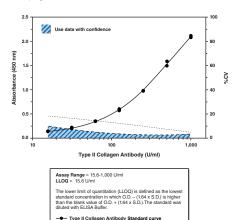
#### Monoacylglycerol Lipase Inhibitor Screening Assay Kit

- Screen for inhibitors of monoacylglycerol lipase (MAGL)
- Assay 45 samples in duplicate
- Includes JZL 195 as a positive control
- Plate-based, colorimetric measurement (405-415 nm)



#### Mouse Anti-Type II Collagen IgG Assay Kit (bovine) 500410

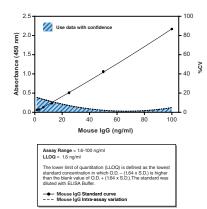
- Measure mouse anti-bovine type II collagen (anti-CII) IgG in mouse plasma or serum
- Employ an affinity-purified anti-CII polyclonal antibody standard isolated from mice with collagen-induced arthritis
- Assay Range: 15.6-1,000 U/ml with an LLOQ of 15.6 U/ml
- Rapid assay; get results in under 4 hours



#### Mouse IgG ELISA Kit

501240

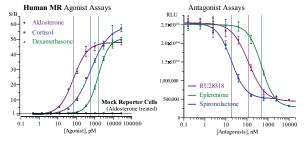
- Measure total mouse IgG from plasma, serum, ascites, and cell culture supernatant
- Convenient 96-well plate format
- Assay Range: 1.6-100 ng/ml with an LLOQ of 1.6 ng/ml
- Get accurate results in under 3 hours



#### MR (human) Reporter Assay Kit

15740

- A nuclear receptor cell-based reporter assay
- · See www.caymanchem.com for more details

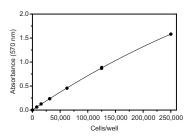




#### MTT Cell Proliferation Assay Kit

10009365

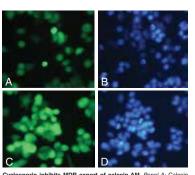
- A colorimetric, non-radioactive assay for assessing cell viability and proliferation
- MTT is reduced to insoluble, purple formazan by NAD(P)Hoxidoreductases in living cells
- Detect 10,000-250,000 cells per well



#### Multi-Drug Resistance Assay Kit (Calcein AM)

600370

- A non-radioactive assay for multi-drug resistance associated protein (MRP) function in living cells
- Employ calcein AM, a substrate for multidrug resistance proteins, including P-gp and MRP
- Cyclosporin A and verapamil are included as positive controls of MDR inhibition
- Assess by flow cytometry, fluorescence microscopy, or plate reader

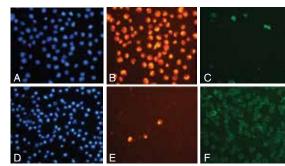


Hoechst Dve staining of MES-SA/Dx5 cells co Panel A. Panel C: Calcein AM staining of MES-SA/Dx5 cells treated with 20 µM Cyclosporin A. Panel D: Hoechst Dye staining of MES-SA/Dx5 cells corresponding to those in Panel C.

#### Multi-Parameter Apoptosis Assay Kit

600330

- Four different reagents for monitoring distinct events in apoptosis
- Contains Annexin V, TMRE, 7-AAD, and Hoechst Dye
- Sufficient reagents to analyze 100 samples for flow cytometry or 500 samples in a 96-well plate format
- 1 hour staining protocol

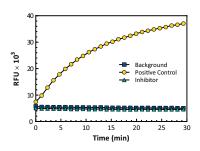


Staurosporine induced apoptosis in Jurkat cells. Jurkat cells treated with vehicle (control) or 2.5 µg/ml staurosporine (treatment) were stained with Hoechst Dye, TMRE or Annexin V FITC. Panels A-C: Control group. Panels D-F: Staurosporine-treated group. Panels A and D: Hoechst staining. Panels B and E: TMRE staining. Panels C and F: Annexin V FITC staining.

#### Myeloperoxidase Chlorination Fluorometric Assay Kit

10006438

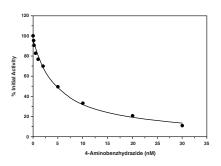
- Measure myeloperoxidase (MPO) chlorination activity in cell lysates and purified preparations
- Assay 37 samples in duplicate
- Plate-based fluorometric measurement (ex 480-495 nm, em 515-525 nm)



#### Myeloperoxidase Inhibitor Screening Assay Kit

700170

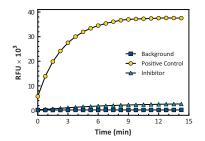
- Screen for inhibitors of myeloperoxidase (MPO)
- Includes reagents to measure both the chlorination and peroxidation activities
- Assay 45 samples in duplicate
- Plate-based, fluorometric measurement (ex 480-495 nm, em 515-525 nm, ex 530-540 nm, em 585-595 nm)



#### Myeloperoxidase Peroxidation Fluorometric Assay Kit

700160

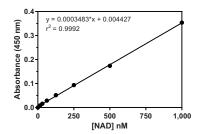
- Measure myeloperoxidase (MPO) peroxidase activity in cell lysates and purified preparations
- Assay 37 samples in duplicate
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### NAD/NADH Cell-Based Assay Kit

• Utilize cellular nicotinamide adenine dinucleotide (NAD+) as a cofactor in a chromogenic enzymatic reaction

- A NAD+ standard curve allows for accurate quantification of cellular NAD levels
- Rapid assay; get results in under 2 hours

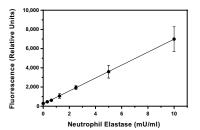


#### Neutrophil Elastase Activity Assay Kit

600610

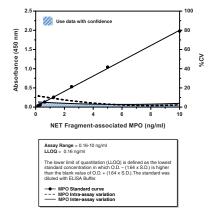
600480

- Neutrophils isolated from whole blood (protocol included)
- Measure neutrophil elastase released by neutrophils after various experimental conditions
- Employ the specific elastase substrate (Z-Ala-Ala-Ala-Ala)2Rh110
- Includes reagents to isolate neutrophils from whole blood and PMA to stimulate elastase release
- Assay performed in provided black 96-well plates
- Neutrophil (mouse) Isolation Kit (Item No. 601070) available to isolate mouse neutrophils from peritoneal lavage or bone marrow



#### Neutrophil Extracellular Trap (MPO-Histone) ELISA Kit

- Detect fragments of neutrophil extracellular traps (NETs) that remain after incomplete digestion and clearance from human samples
- A sandwich ELISA using antibodies specific for MPO and citrullinated histone H3
- Quantitative; results are expressed in ng/ml of NET-associated MPO



Neutrophil Extracellular Trap (NET) Assay Kit

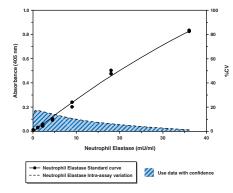
NF-κB (human p50) Transcription 601010

Factor Assay Kit

10006912

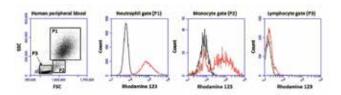
• Induce and detect NET formation in vivo

- Elastase-based readout
- Non-dsDNA readout eliminates false positives from DNA
- Adaptable to multiple species
- Neutrophil (mouse) Isolation Kit (Item No. 601070) available to isolate mouse neutrophils from peritoneal lavage or bone marrow



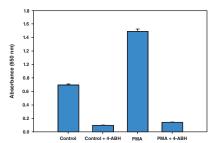
#### Neutrophil/Monocyte Respiratory Burst Assay Kit 601130

- · Samples include whole blood from any species or cells of any type that are capable of producing a nicotinamide adenine dinucleotide, reduced (NADPH) oxidase-dependent respiratory burst
- Induce and quantify a respiratory burst response
- Analyze by flow cytometry



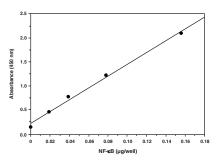
#### Neutrophil Myeloperoxidase Activity Assay Kit 600620

- Assay the release of enzymatically-active myeloperoxidase (MPO) from activated phagocytes
- Utilize TMB as a chromogenic substrate for MPO
- Includes a specific inhibitor of MPO function to verify specificity
- Includes reagents needed to isolate neutrophils from human whole



# • A sensitive, non-radioactive method of detecting NF-kB (p50)

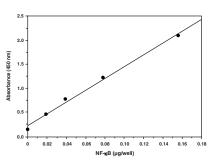
- from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



#### NF-κB (p65) Transcription Factor Assay Kit

10007889

- A sensitive, non-radioactive method of detecting NF-kB (p65) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



#### Nitrate/Nitrite Colorimetric Assay Kit

Nitrate/Nitrite Colorimetric Assay Kit (LDH method)

> serum, and tissue homogenates • Assay 41 samples in duplicate

• Assay Range: 5-25 μM (nitrate and/or nitrite) • Plate-based, colorimetric measurement (530-550 nm)

• Measure nitrate, nitrite, and/or total nitrate/nitrites in

plasma/serum, urine, tissue culture media, and tissue homogenates

• Assay 26 samples in triplicate or 40 samples in duplicate/96-well plate

• Designed to measure nitric oxide synthase (NOS) activity in vitro • Can also be used to measure nitrate and nitrite in urine, plasma,

#### NOS Activity Assay Kit

780001

760871

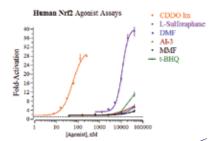
#### 781001

- Measure nitric oxide synthase (NOS) activity in cell lysates and purified preparations
- Material for 50 reactions
- Radioactive substrate not supplied

#### Nrf2 (human) Reporter Assay Kit

17734

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details





#### Nrf2 Transcription Factor Assay Kit

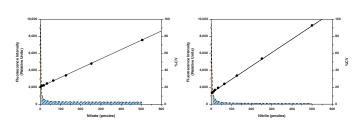
600590

- A sensitive, non-radioactive method of detecting nuclear factor E2related factor 2 (Nrf2) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays (EMSAs)
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates

# Nitrate/Nitrite Fluorometric Assay Kit

780051

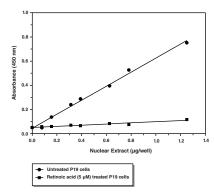
- Measure nitrate, nitrite, and/or total nitrate/nitrites in plasma, serum, tissue culture media, and tissue homogenates
- Assay 40 samples in duplicate/96-well plate
- Assay Range: 60 nM-3.85 μM (nitrate and/or nitrite)
- Plate-based, fluorometric measurement (ex 360-365 nm, em 430 nm)



#### S-Nitrosylated Protein Detection Kit

- A modified 'Biotin-switch' assay
- Direct visualization of S-nitrosylated (S-NO) proteins in whole cells or tissues
- Colorimetric or fluorometric detection

- A sensitive, non-radioactive method of detecting Oct4 from whole
- 96-well plate format replaces electrophoretic mobility shift assays (FMSAs)
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates

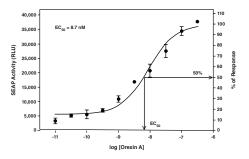


#### Orexin 1 Receptor Reporter Assay Kit

600240

601080

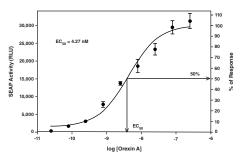
- Reverse transfection reporter assay to screen for orexin 1 receptor (OX1R) agonists
- Assay 46 samples in duplicate
- Includes orexin A as a positive control
- · Plate-based, luminometric measurement



#### Orexin 2 Receptor Reporter Assay Kit

#### 600250

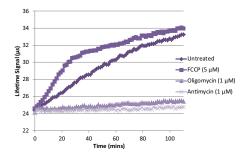
- A 96-well plate coated with DNA encoding both orexin 2 receptor (OX2R) and a cAMP response element-regulated SEAP
- Cells grown on the complex will express both OX2R and the SEAP
- Binding of OX2R agonists induces expression of SEAP
- SEAP secreted into the supernatant cleaves a luminescence-based substrate (included)



#### Oxygen Consumption Rate Assay Kit (MitoXpress®-Xtra HS Method)

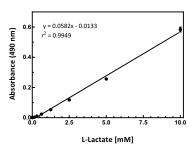
600800

- Measure oxygen consumption rate (OCR) without the need for an oxygen electrode
- Utilize MitoXpress®-Xtra, a phosphorescent oxygen probe
- Includes antimycin A, an inhibitor of oxygen consumption, as a control
- Includes glucose oxidase as a reference for oxygen depletion



#### Oxygen Consumption/Glycolysis Dual Assay Kit

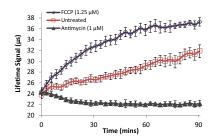
- Measure both oxygen consumption and glycolysis
- Utilize MitoXpress®-Xtra, a phosphorescent oxygen probe
- Includes antimycin A, an inhibitor of oxygen consumption, as a control



#### Oxygen Consumption/MitoMembrane Potential Dual Assay Kit

600880

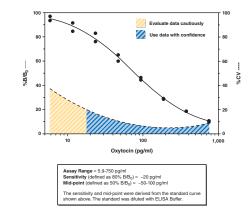
- Measure both oxygen consumption rate (OCR) and mitochondrial membrane potential
- Utilize MitoXpress®-Xtra, a phosphorescent oxygen probe
- Employ the cationic dye, JC-1, to determine mitochondrial membrane potential
- Includes antimycin A, an inhibitor of oxygen consumption, as a control
- Includes glucose oxidase as a reference for oxygen depletion



#### • Measure oxytocin in plasma

Oxytocin ELISA Kit

- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure oxytocin levels down to 20 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

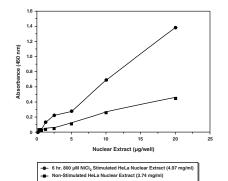


#### p53 Transcription Factor Assay Kit

600020

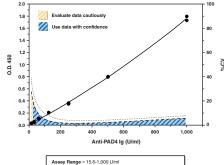
500440

- A sensitive, non-radioactive method of detecting p53 from whole
- 96-well plate format replaces electrophoretic mobility shift assays
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



#### • Measure anti-protein arginine deiminase 4 (anti-PAD4)

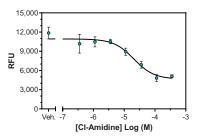
- autoantibodies of any isotype (IgM, IgG, IgA) • For use in human plasma and serum (no need to purify samples)
- Assay Range: 15.6-1,000 U/ml with an LLOQ of 31.3 U/ml





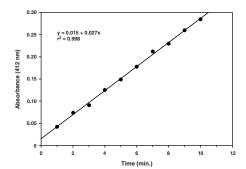
#### PAD4 Inhibitor Screening Assay Kit (Ammonia)

- Screen for inhibitors of protein arginine deiminase 4 (PAD4)
- Assay 45 samples in duplicate
- Plate-based, fluorometric measurement (ex 405-415 nm, em 470-480 nm)



#### PAF Acetylhydrolase Assay Kit

- Measure PAF acetylhydrolase (PAF-AH) activity in plasma, serum, tissue homogenates, and cell lysates
- · Assay 23 samples in duplicate
- Measure PAF-AH activity down to 0.02 U/ml
- Plate-based, colorimetric measurement (405-415 nm)

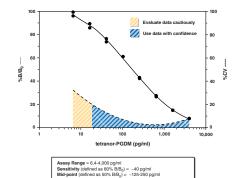


- Screen for inhibitors of PAF acetylhydrolase (PAF-AH)
- Assay 45 samples in duplicate
- Plate-based, colorimetric measurement (405-415 nm)

#### tetranor-PGDM ELISA Kit

#### 501001

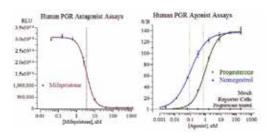
- Detect tetranor-prostaglandin D metabolite (tetranor-PDGM), a major metabolite of prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) in urine
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure tetranor-PGDM levels down to 40 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



#### PGR (human) Reporter Assay Kit

15759

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

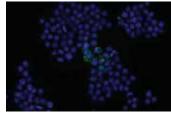




#### Phagocytosis Assay Kit (IgG FITC)

500290

- Detect phagocytosis of rabbit IgG-coated latex beads
- Validated in human peripheral blood monocytes and mouse peritoneal macrophages
- Phagocytosis detectable in 1-4 hours at 37°C
- Assess by flow cytometry or fluorescence microscopy

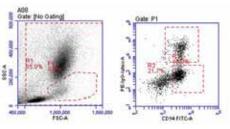


RAW 264.7 cells phagocytose opsonized particles.

- Detect phagocytosis of rabbit IgG-coated latex beads
- Validated in human peripheral blood monocytes and mouse peritoneal macrophages
- Phagocytosis detectable in 1-4 hours at 37°C

Phagocytosis Assay Kit (IgG PE)

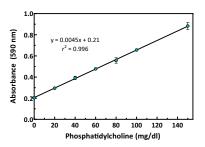
Assess by flow cytometry or fluorescence microscopy



Peripheral blood monocytes (CD14+) phagocytose opsonized particles

#### Phosphatidylcholine Colorimetric Assay Kit 10009926

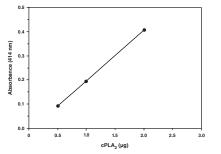
- Measure phosphatidylcholine (PC) in serum or plasma
- · Assay 41 samples in duplicate
- Assay Range: 20-150 mg/dl
- Plate-based, colorimetric measurement (580-600 nm)



#### cPLA<sub>2</sub> Assay Kit

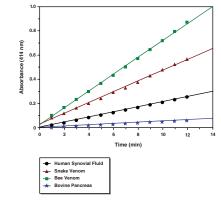
765021

- Measure calcium-dependent cytosolic phospholipase A<sub>2</sub> (cPLA<sub>2</sub>) activity in cell lysates and tissue homogenates that are known to only contain cPLA<sub>2</sub>, or in purified enzyme preparations
- Assay 46 samples in duplicate
- Measure cPLA<sub>2</sub> activity down to 0.0035 U/ml
- Plate-based, colorimetric measurement (405-420 nm)



#### • Measure secretory phospholipase A2 (sPLA2) activity

- · Assay 46 samples in duplicate
- Measure sPLA, activity down to 0.02 U/ml
- Plate-based, colorimetric measurement (405-414 nm)

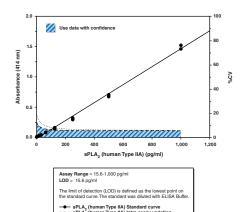


#### sPLA<sub>2</sub> (human Type IIA) ELISA Kit

585000

765001

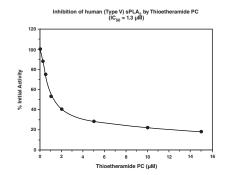
- Measure secretory phospholipase A2 (sPLA2) levels in plasma, synovial fluid, and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Assay Range: 15.6-1,000 pg/ml with an LOD of 15.6 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



#### sPLA<sub>2</sub> (Type V) Inhibitor Screening Assay Kit

10004883

- Screen for inhibitors of Type V secretory phospholipase A2 (sPLA2)
- Assay 45 samples in duplicate
- Plate-based, colorimetric measurement (405-420 nm)



#### PPAR (human) Reporter Assays Panel

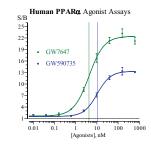
- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



#### PPARα (human) Reporter Assay Kit

15730

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

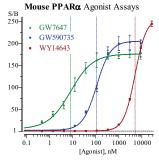




#### PPARα (mouse) Reporter Assay Kit

15762

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



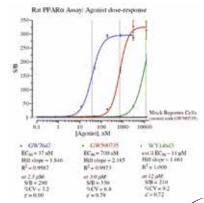


15764

#### PPARα (rat) Reporter Assay Kit

A nuclear receptor cell-based reporter assay

• See www.caymanchem.com for more details





PPARα Transcription Factor Assay Kit

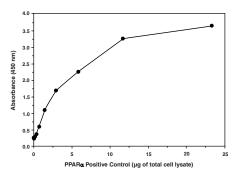
10006915

PPARδ (human) Reporter Assay Kit

PPARδ Transcription Factor Assay Kit

PPARy Transcription Factor Assay Kit

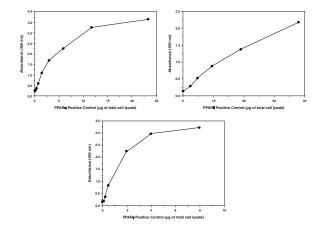
- A sensitive, non-radioactive method of detecting peroxisome proliferator-activated receptor α (PPARα) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays (FMSAs)
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



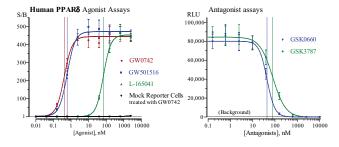
#### PPAR $\alpha$ , $\delta$ , $\gamma$ Complete Transcription Factor Assay Kit

10008878

- A sensitive, non-radioactive method of detecting peroxisome proliferator-activated receptor (PPAR)α, PPARγ, and PPARδ/β from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Single 96-well plate that measures all three PPAR isoforms
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



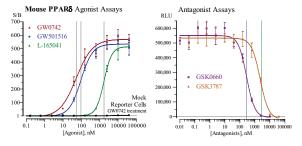


#### PPARδ (mouse) Reporter Assay Kit

15763

15731

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

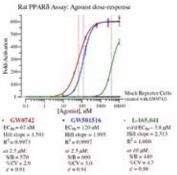




#### PPARδ (rat) Reporter Assay Kit

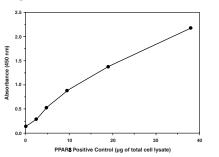
15765

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details





- A sensitive, non-radioactive method of detecting peroxisome proliferator-activated receptor  $\delta$  (PPAR $\delta$ ) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays (EMSAs)
- Capture the transcription factor using a specific dsDNA sequence
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Plate-based, colorimetric measurement (405-420 nm)
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates

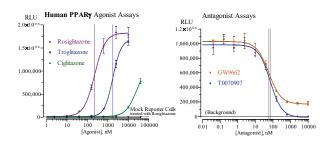


#### PPARγ (human) Reporter Assay Kit

15729

10006914

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

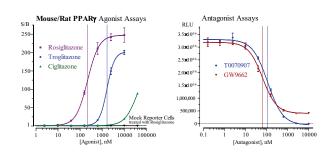




#### PPARy (mouse/rat) Reporter Assay Kit

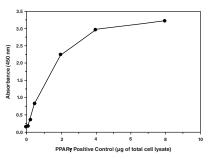
15761

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details





- A sensitive, non-radioactive method of detecting peroxisome
- proliferator-activated receptor γ (PPARγ) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays (EMSAs)
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



#### Pregnane X (human) Receptor Assay Kit

17528

17585

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



#### Pregnane X (rat) Receptor Assay Kit

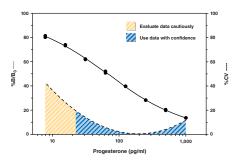
• A nuclear receptor cell-based reporter assay

• See www.caymanchem.com for more details



#### Progesterone ELISA Kit

- Measure progesterone in plasma and other sample matrices
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure progesterone levels down to 10 pg/ml
- Rapid assay; get results in under 4 hours



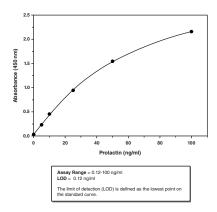
Prolactin (human) ELISA Kit

#### Prostaglandin D<sub>2</sub> ELISA Kit

#### Prostaglandin D<sub>2</sub>-MOX ELISA Kit

#### Prostaglandin E Metabolite ELISA Kit

- Measure prolactin in human serum and plasma
- Assay 27 samples in triplicate or 41 samples in duplicate
- Assay Range: 0.12-100 ng/ml with an LOD of 0.12 ng/ml
- Rapid assay; get results in under 2 hours

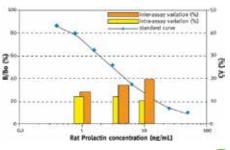


#### Prolactin (rat) EIA Kit

589701

500730

- · Measure prolactin in rat plasma, serum, and cell culture supernatant samples
- Assay 33 samples in duplicate
- Incubation: Overnight
- No extraction necessary

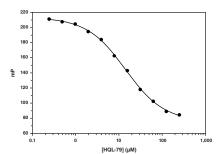


# SPI DIO DELLO

#### Prostaglandin D Synthase (hematopoietic-type) FP-Based Inhibitor Screening Assay Kit - Green

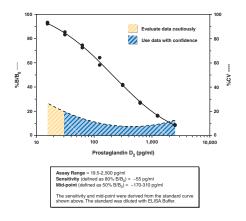
600007

- Screen for inhibitors of prostaglandin D synthase (hematopoietic-type) (H-PGDS)
- Designed for high-throughput screening
- Includes HQL-79 as a positive control
- Plate-based, fluorescence polarization measurement (ex 470 nm, em 530 nm)



#### • Measure prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) in multiple sample types

- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure PGD, levels down to 55 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

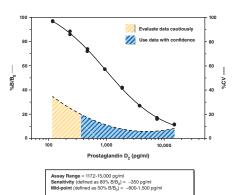


#### Prostaglandin D<sub>2</sub> Express ELISA Kit

512041

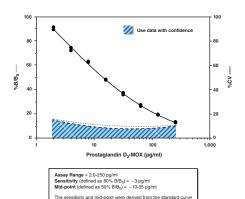
512031

- Measure prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) in multiple sample types
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure PGD, levels down to 350 pg/ml
- Rapid assay; get results in under 4 hours



#### • Treatment of samples with methoxylamine (MOX) hydrochloride converts unstable prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) to stable PGD<sub>2</sub>-MOX

- Highly sensitive
- Measure PGD<sub>2</sub>-MOX levels down to 3 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

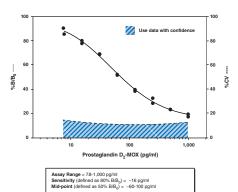


#### Prostaglandin D<sub>2</sub>-MOX Express ELISA Kit

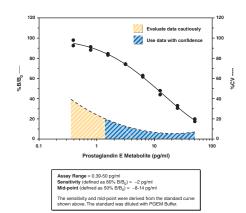
#### 500151

512011

- Treatment of samples with methoxylamine (MOX) hydrochloride converts unstable prostaglandin D<sub>2</sub> (PGD<sub>2</sub>) to stable PGD<sub>2</sub>-MOX; MOX derivatization reagent included with assay
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure PGD<sub>2</sub>-MOX levels down to 16 pg/ml
- Rapid assay; get results in under 4 hours

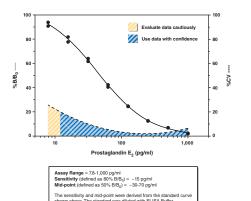


- Converts major urinary and plasma metabolites of prostaglandin E<sub>a</sub> (PGE<sub>2</sub>) to a single, stable derivative that is easily quantified
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure PGEM levels down to 2 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm

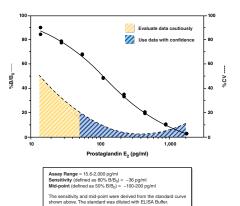


#### Prostaglandin E2 ELISA Kit - Monoclonal

- Measure prostaglandin E<sub>2</sub> (PGE<sub>2</sub>) in multiple sample types
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure PGE, levels down to 15 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



- Measure prostaglandin E<sub>2</sub> (PGE<sub>2</sub>) in multiple sample types
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure PGE, levels down to 36 pg/ml
- Rapid assay; get results in under 4 hours

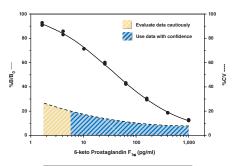


#### 6-keto Prostaglandin F<sub>10</sub> ELISA Kit

515211

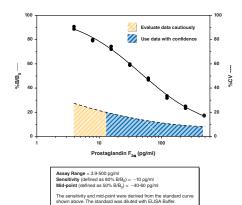
500141

- Detect 6-keto prostaglandin F<sub>1a</sub> (PGF<sub>1a</sub>), a surrogate for prostacyclin synthesis in multiple sample types
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 6-keto PGF<sub>1g</sub> levels down to 6 pg/ml
- Incubation: 18 hours Development: 90-120 minutes Read: Colorimetric at 405-420 nm



Assay Range = 1.6-1,000 pg/ml Sensitivity (defined as 80% B/B<sub>0</sub>) = -6 pg/ml Mid-point (defined as 50% B/B<sub>0</sub>) = -20-60 pg/ml sensitivity and mid-point were derived from the standar in above. The standard was diluted with ELISA Buffer.

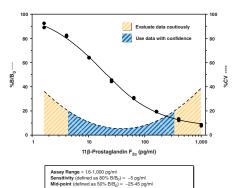
- Measure prostaglandin  $F_{2a}$  (PGF $_{2a}$ ) in multiple sample types
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure PGF<sub>20</sub> levels down to 10 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



#### 11β-Prostaglandin F<sub>2a</sub> ELISA Kit

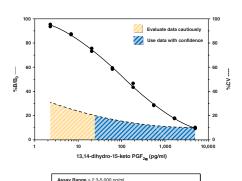
516521

- Detect 11β-prostaglandin F<sub>20</sub> (PGF<sub>20</sub>), the primary metabolite of PGD<sub>2</sub>, in plasma or urine
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 11β-PGF<sub>2a</sub> levels down to 5 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



#### 13,14-dihydro-15-keto Prostaglandin F<sub>20</sub> ELISA Kit 516671

- Detect 13,14-dihydro-15-keto prostaglandin F<sub>20</sub> (PGF<sub>20</sub>), a stable metabolite of PGF
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 13,14-dihydro-15-keto PGF<sub>20</sub> levels down to 15 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

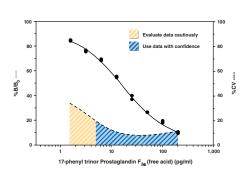


### Assay Range = 2.3-5,000 pg/ml Sensitivity (defined as $80\% B/B_0$ ) = -15 pg/mlMid-point (defined as $50\% B/B_0$ ) = -90-150 pg/mThe sensitivity and mid-point were derived from the standard shown above. The standard was diluted with ELISA Buffer.

#### 17-phenyl trinor Prostaglandin F<sub>2a</sub> ELISA Kit

516821

- Measure 17-phenyl trinor prostaglandin F<sub>2</sub> (PGF<sub>2</sub>) (also known as bimatoprost)
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 17-phenyl trinor PGF<sub>20</sub> levels down to 2.6 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



Assay Range = 1.56-200 pg/ml Sensitivity (defined as 80% B/B<sub>0</sub>) = -2.6 pg/ml Mid-point (defined as 50% B/B<sub>0</sub>) = -15 pg/ml

# 1.000

• Measure prostaglandin I metabolites (PGIMs) in urine

• Measure PGIM down to 120 pg/ml

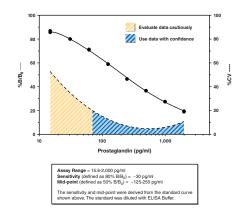
Convenient overnight format

Assay Range = 39-5,000 pg/ml Sensitivity (defined as 80% B/B<sub>0</sub>) =  $\sim$ 120 pg/ml Mid-point (defined as 50% B/B<sub>0</sub>) =  $\sim$ 500 pg/ml The sensitivity and mid-point were derived from the standar shown above. The standard was diluted with Assay Buffer.

#### Prostaglandin Screening ELISA Kit

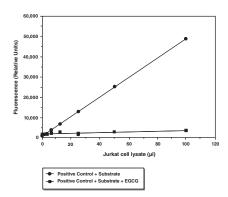
514012

- Measure total prostaglandin (PG) levels in tissue culture medium
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure total PG levels down to 30 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



#### 20S Proteasome Assay Kit

- A functional assay for detecting activity of the 20S proteasome in
- Utilize SUC-LLVY-AMC as a specific substrate for the 20S proteasome
- Includes a Jurkat cell lysate as a positive control
- Includes EGCG, a specific 20S proteasome inhibitor



Protein Carbonyl Colorimetric Assay Kit

Rabbit Anti-Goat Fluorophore Selection Kit

16544

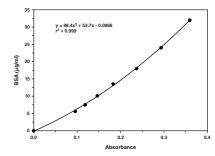
- Measure protein carbonyl content in plasma, serum, urine, tissue homogenates, and cell lysates
- Assay 24 samples in duplicate
- Plate-based, colorimetric measurement (360-385 nm)

#### Protein Determination Kit

704002

10005020

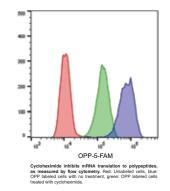
- Measure total protein concentration
- Assay Range: 5.6-32 μg/ml
- Plate-based, colorimetric measurement (595 nm)



#### Protein Synthesis Assay Kit

601100

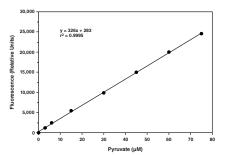
- Non-radioactive, simple procedure
- Measure nascent protein synthesis in cells
- O-Propargyl-puromycin (OPP) and 5 FAM-Azide used to detect nascent proteins by flow cytometry, microscopy, or plate reader
- Sufficient reagents to run up to 100 samples



#### Pyruvate Assay Kit

700470

- Measure pyruvate in serum, plasma, blood, urine, and saliva
- Assay 40 samples in duplicate
- Assay Range: 3-70 μM
- Plate-based, fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### • Comprehensive: Columbia Biosciences' fluorophore selection kits include the secondary antibody individually conjugated to seven unique fluorophores

- Efficient: The ability to evaluate the various labeled conjugates side-by-side allows scientists to rapidly optimize their reagents
- Cost-effective: Order one kit and get seven conjugated secondary antibodies. The fluorophore selection kit reduces costs associated with purchasing and shipping full-size antibody conjugates.





#### RAR (human) Reporter Assays Panel

15755

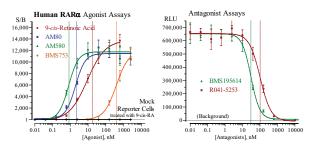
- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



#### RARα (human) Reporter Assay Kit

15754

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

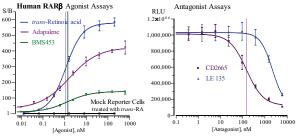




#### • A nuclear receptor cell-based reporter assay

RARβ (human) Reporter Assay Kit

• See www.caymanchem.com for more details

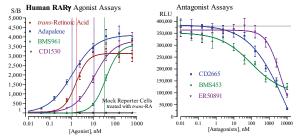




#### RARy (human) Reporter Assay Kit

#### 15752

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

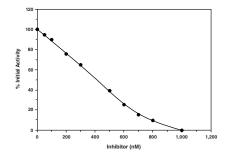




#### Renin Inhibitor Screening Assay Kit

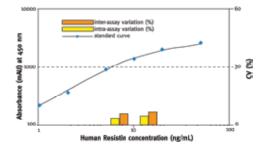
#### 10006270

- Screen for inhibitors of renin aspartyl protease activity
- Assay 45 samples in duplicate
- Plate-based, fluorometric measurement (ex 335-345 nm, em 484-510 nm)



#### Resistin (human) EIA Kit

- Measure resistin in human serum, plasma, and culture supernatant
- Assay 41 samples in duplicate
- Incubation: 3 x 1 hour

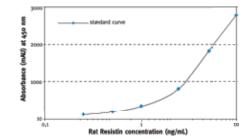




#### Resistin (rat) EIA Kit

10007612

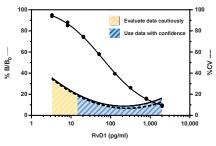
- Measure resistin in mouse serum, plasma, and culture supernatant
- Assay 40 samples in duplicate
- Incubations: 2 x 1 hour | 30 minutes

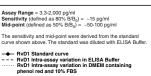




#### Resolvin D1 ELISA Kit

- Detect resolvin D1 (RvD1) in multiple sample types
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure RvD1 levels down to 15 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm





Resolvin D2 ELISA Kit

• Detect resolvin D2 (RvD2) in multiple sample types

• Measure RvD2 levels down to 10 pg/ml

Read: Colorimetric at 405-420 nm

• Assay 24 samples in triplicate or 36 samples in duplicate

• Incubation: 18 hours | Development: 90-120 minutes |

Assay Range = 1.6-1,000 pg/ml Sensitivity (defined as 80% B/B $_0$ ) = -10 pg/ml Mid-point (defined as 50% B/B $_0$ ) = -20-60 pg/ml

Human RORy Inverse-Agonist Assays

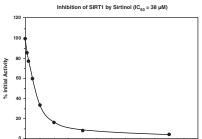
100 Resolvin D2 (pg/ml)

#### RXRα (human) Reporter Assay Kit

#### 15743

#### SIRT1 Direct Fluorescent Screening Assay Kit 10010401

- Assay 45 samples in duplicate
- Plate-based, fluorometric measurement (ex 340-360 nm, em 450-465 nm)

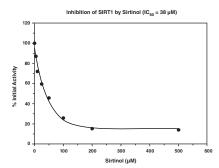


#### SIRT1 FRET-Based Screening Assay Kit

10010991

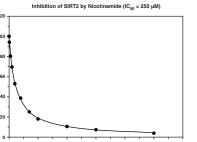
700280

- Screen for modulators of SIRT1 activity
- Assay 45 samples in duplicate
- em 440-465 nm)

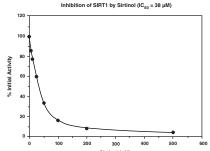


#### SIRT2 Direct Fluorescent Screening Assay Kit

- Screen for modulators of SIRT2 activity
- Assay 45 samples in duplicate
- em 450-465 nm)

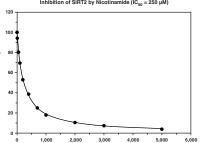


- Screen for modulators of SIRT1 activity



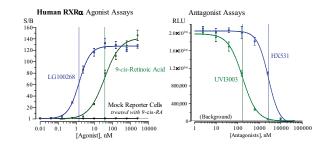
- Plate-based, fluorometric measurement (ex 335-345 nm,

- Plate-based, fluorometric measurement (ex 340-360 nm,



501120

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details





#### RXRβ (human) Reporter Assay Kit

Human RXRB Agonist Assays

• A nuclear receptor cell-based reporter assay

• See www.caymanchem.com for more details

15744

- Semi-Quantitative Biomarker
  - **ELISA Component Kit (anti-mouse)**

Secreted Alkaline Phosphatase

Reporter Gene Assay Kit (Luminescence)

• Rapid assay; get results within 1 hour

phosphatase (SEAP) from transfected cells

• A simple chemiluminescence assay for measuring secreted alkaline

• Detect SEAP activity in culture medium in the mU/well range

• Contains enough reagents to run three 96-well plates

- A generic assay for measurement of fish biomarkers
- See www.caymanchem.com for more details

bisense

10008659

10007680

600260

#### Semi-Quantitative Biomarker ELISA Component Kit (anti-rabbit)

- A generic assay for measurement of fish biomarkers
- See www.caymanchem.com for more details

bisense

600490

# **MINDIGO**

# **MINDIGO**

#### RXR (human) Reporter Assays Panel

RORγ (human) Reporter Assay Kit

• A nuclear receptor cell-based reporter assay

• See www.caymanchem.com for more details

#### 15746

15757

• A nuclear receptor cell-based reporter assay

- TO901317

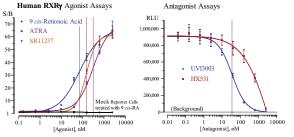
SR 1001

• See www.caymanchem.com for more details



#### RXRγ (human) Reporter Assay Kit

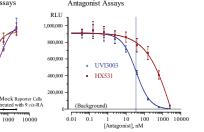
1.2×10





#### SET7/9 SAM-Screener™ Assay Kit 15745

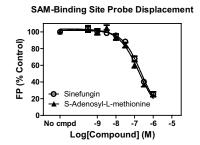
- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details



Antagonist Assays

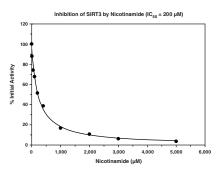


- Homogeneous, fluorescence polarization assay to screen for SAMbinding site inhibitors
- Designed for high-throughput screening
- Plate-based, fluorescence polarization measurement (ex 575 nm, em 620 nm)



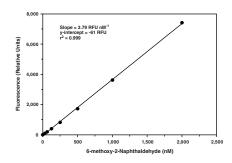
SIRT3 Direct Fluorescent Screening Assay Kit 10011566

- · Screen for modulators of SIRT3 activity • Assay 45 samples in duplicate
- Plate-based, fluorometric measurement (ex 340-360 nm, em 450-465 nm)



#### Soluble Epoxide Hydrolase Cell-Based Assay Kit 600090

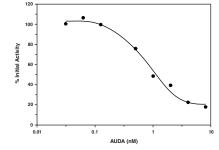
- A convenient 96-well fluorecence-based assay for detecting soluble epoxide hydrolase (sEH) activity in whole cells
- Utilize epoxy fluor 7, a sensitive fluorescent substrate for sEH
- Quantifiable using the included 6-methoxy-2-naphthaldehyde standard
- Includes AUDA, a selective epoxide hydrolase inhibitor, for checking specificity of the reaction



#### Soluble Epoxide Hydrolase Inhibitor Screening Assay Kit

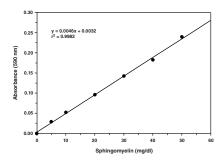
10011671

- Screen for inhibitors of soluble epoxide hydrolase (sEH)
- Assay 45 samples in duplicate
- Plate-based, fluorometric measurement (ex 330 nm, em 465 nm)



### Sphingomyelin Colorimetric Assay Kit

- Measure sphingomyelin (SM) in plasma and serum
- Assay 41 samples in duplicate • Assay Range: 5-50 mg/dl
- Plate-based, colorimetric measurement (585-600 nm)

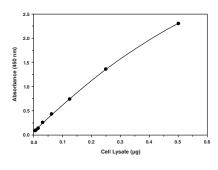


#### SREBP-1 Transcription Factor Assay Kit

10010854

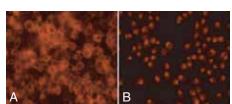
10009928

- A sensitive, non-radioactive method of detecting sterol regulatory element-binding protein-1 (SREBP-1) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays (EMSAs)
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates



#### SREBP-2 Cell-Based Translocation Assay Kit 10009239

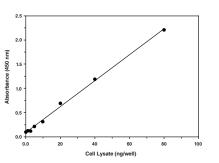
- Detect translocation of sterol regulatory element-binding protein-2 (SREBP-2) into the nucleus of treated and fixed cells by immunofluorescence
- Employ a highly specific antibody to SREBP-2
- An inhibitor of cholesterol trafficking is included as a positive control



Translocation of SREBP-2 into nuclei by 24 µM U-18666A.

#### SREBP-2 Transcription Factor Assay Kit

- A sensitive, non-radioactive method of detecting sterol regulatory element-binding protein-2 (SREBP-2) from whole cell lysates
- 96-well plate format replaces electrophoretic mobility shift assays (FMSAs)
- Capture the transcription factor using a specific dsDNA sequence bound to the plate
- Detect the dsDNA-bound transcription factor with specific antibodies in an ELISA format
- Nuclear Extraction Kit (Item No. 10009277) available to aid in the isolation of nuclear and cytoplasmic fractions from cell lysates or tissue homogenates

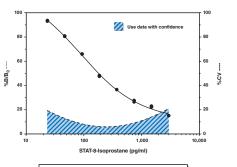


#### STAT-8-Isoprostane ELISA Kit

500431

10007819

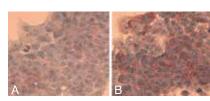
- Measure 8-isoprostane, a biomarker of oxidative stress
- Assav 24 samples in triplicate or 36 samples in duplicate
- Measure 8-isoprostane levels down to 45 pg/ml
- Extremely rapid assay; get results in about 2 hours
- Alkaline phosphatase used as tracer



#### Steatosis Colorimetric Assay Kit

10012643

- Detect excessive lipid accumulation in cells
- Assess by microscopy or plate reader
- An inducer of lipid accumulation is included as a positive control

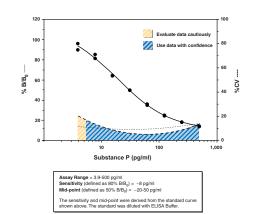


Effect of chloroquine on lipid droplet accumulation in HepG2

#### Substance P ELISA Kit

• Measure substance P in multiple sample matricies from multiple

- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure substance P levels down to 8 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm

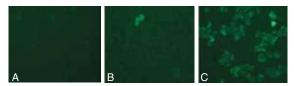


#### Sulfenylated Protein Cell-Based Detection Kit

600320

583751

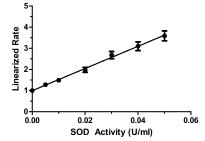
- Contains the cell-permeable probe DAz-2 to detect sulfenic acidmodified proteins in cells
- Includes epigallocatechin gallate (EGCG) as a positive control
- Detect by flow cytometry, fluorescence microscopy, or plate reader



Hydrogen peroxide induces protein sulfenylation in RAW 264.7 cells. Panel A: Control, no DAz-2. Panel B: Cells treated with vehicle. Panel C: Cells treated with 0.25% H.O.

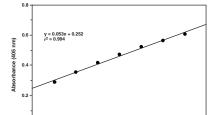
#### Superoxide Dismutase Assay Kit

- Measure copper/zinc, manganese, and iron superoxide dismutase (SOD) in tissues, plasma, serum, and cell lysates
- Assay 41 samples in duplicate
- Measure SOD activity down to 0.005 U/ml



• Measure thioredoxin reductase (TrxR) activity in tissue homogenates

- - and cell lysates • Assay 22 samples in duplicate
  - Measure TrxR down to 0.08 U/ml
  - Plate-based colorimetric measurement (405-414 nm)



#### Thioredoxin Reductase Fluorescent Activity Assay Kit

11529

- Convenient 96-well plate format
- Measure thioredoxin reductase (TrxR) in cell lysates and tissue homogenates
- Active TrxR and human Trx-1 provided with kit



11526

#### Thioredoxin/Thioredoxin Reductase Mammalian Assay Kit

- Convenient 96-well plate format
- Measure both thioredoxin reductase (TrxR) and Trx in cell lysates and tissue homogenates
- Active TrxR and human Trx-1 provided with kit

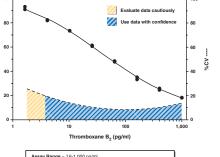
#### IMCO

501020

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

• Measure thromboxane B<sub>2</sub> (TXB<sub>2</sub>) in multiple sample types

- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure TXB, levels down to 5 pg/ml
- Incubation: 18 hours | Development: 90-120 minutes | Read: Colorimetric at 405-420 nm



• A complete kit for conjugating APC to IgG





#### SureLight® APC Conjugation Kit 2 mg IgG

- A complete kit for conjugating APC to IgG
- See www.caymanchem.com for more details



#### SureLight® R-PE Conjugation Kit

13436

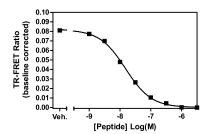
- A complete kit for conjugating R-PE to proteins
- See www.caymanchem.com for more details



#### TAF1 bromodomain 1 TR-FRET Assay Kit

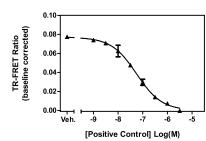
600870

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- · Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### TAF1 bromodomains 1 and 2 TR-FRET Assay Kit 600930

- Homogeneous, time-resolved fluorescence energy transfer (TR-FRET) assay to screen for inhibitors
- Designed for high-throughput screening
- Plate-based, time-resolved fluorometric measurement (ex 340 nm, em 620 and 670 nm)



#### 13412

13411

• Measure thiobarbituric acid reactive substances (TBARS) in

• Plate-based, colorimetric measurement (530-540 nm) or

fluorometric measurement (ex 530 nm, em 550 nm)

plasma, serum, urine, tissue homogenates, and cell lysates

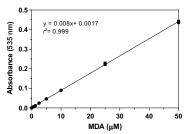


• Assay 40 samples in duplicate

• Assay Range: 0.0625-50 μM

700870

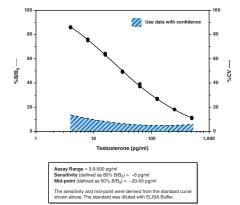
- Measure thiobarbituric acid reactive substances (TBARS) in plasma, serum, urine, tissue homogenates, and cell lysates
- Assay 40 samples in duplicate
- Assay Range: 0.0625-50 μM
- Plate-based, colorimetric measurement (530-540 nm) or fluorometric measurement (ex 530 nm, em 550 nm)



#### Testosterone ELISA Kit

582701

- Measure testosterone in multiple sample types
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure testosterone levels down to 6 pg/ml
- Incubation: 2 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm

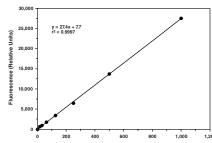


#### Thiol Detection Assay Kit

• Measure free thiol content in plasma, serum, urine, cell lysates,

and tissue homogenates

• Plate-based fluorometric measurement (ex 380-390 nm,



#### Thioredoxin Activity Fluorescent V2 Assay Kit

• Active TrxR and human Trx-1 provided with kit

IMCO



700340

· Assay 40 samples in duplicate

• Assay Range: 15 nM-1 μM

11-keto Testosterone ELISA Kit

• Detect the primary androgenic steroid in fish

Read: Colorimetric at 405-420 nm

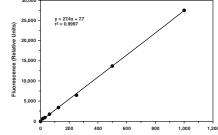
• Assay 24 samples in triplicate or 36 samples in duplicate

• Measure 11-keto testosterone levels down to 1.3 pg/ml

• Incubation: 18 hours | Development: 90-120 minutes |

em 510-520 nm)

• Includes both a glutathione and cysteine standard



# 15493

#### · Convenient 96-well plate format

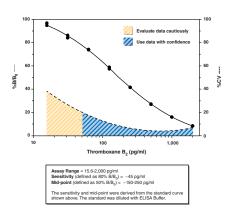
- Measure thioredoxin (Trx) activity in cell lysates and tissue homogenates

10004023

• Measure thromboxane B<sub>2</sub> (TXB<sub>2</sub>) in multiple sample types

• Assay 24 samples in triplicate or 36 samples in duplicate

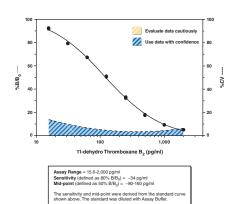
- Measure TXB, levels down to 45 pg/ml
- Rapid assay; get results in under 4 hours



#### 11-dehydro Thromboxane B<sub>2</sub> ELISA Kit - Monoclonal

519510

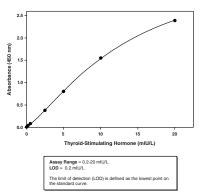
- Detect 11-dehydro thromboxane B2 (TXB2), a stable urinary metabolite of TXA, and TXB,
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure 11-dehydro TXB<sub>2</sub> levels down to 34 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



#### Thyroid-Stimulating Hormone ELISA Kit

500740

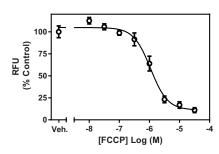
- Measure thyroid-stimulating hormone (TSH) in human serum and plasma
- Assay 26 samples in triplicate or 40 samples in duplicate
- Assay Range: 0.2-20 mlU/L with an LOD of 0.2 mlU/L
- Rapid assay; get results in under 3 hours



#### TMRE Mitochondrial Membrane

Potential Assay Kit 701310

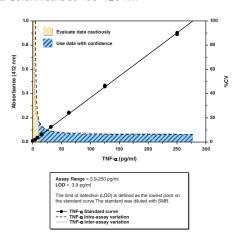
- Detect mitochondrial membrane potential as an indicator of mitochondrial health
- Utilize tetramethylrhodamine ethyl ester (TMRE), a fluorescent, lipophilic, cationic dye
- Staining in less than 1 hour



#### TNF-α (human) ELISA Kit

589201

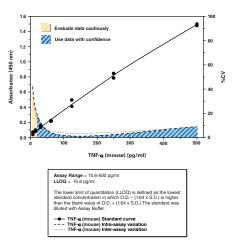
- Measure tumor necrosis factor-α (TNF-α) from human plasma, serum, and other sample matrices
- Assay Range: 3.9-250 pg/ml with an LOD of 3.9 pg/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



#### TNF-α (mouse) ELISA Kit

500850

- Measure tumor necrosis factor-α (TNF-α) from mouse plasma, serum, and other sample matrices
- Assay Range: 15.6-500 pg/ml with an LLOQ of 15.6 pg/ml
- Rapid assay; get results in under 3 hours



#### TR (human) Reporter Assays Panel

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

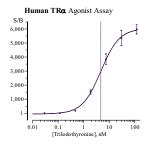


#### TRα (human) Reporter Assay Kit

15749

15751

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

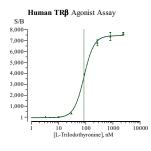


# **MINDIGO**

#### TRβ (human) Reporter Assay Kit

15750

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

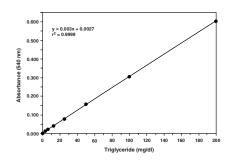




#### Triglyceride Colorimetric Assay Kit

10010303

- Measure triglycerides (TGs) levels in plasma, serum, cell lysates, and tissue homogenates
- Assay 26 samples in triplicate and 40 samples in duplicate
- Assay Range: 3.125 to 200 mg/dl
- Incubation: 15 minutes | Read: Colorimetric at 530-550 nm



#### **Ubiquitin Interact Kit**

• Identify and characterize isolated ubiquitin binding proteins by western blotting or proteomic analysis

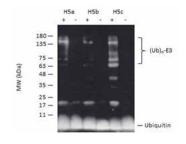
- Demonstrate ubiquitin binding by specific proteins of interest
- Investigate the role of ubiquitin in particular signaling pathways

#### **Ubiquitin Link Kit**

15985

15978

- Develop assays for E3 auto-ubiquitination and substrate ubiquitination
- Determine suitable E2 for use with specific E3 or E3-substrate systems
- Assess E3 ligase activity of novel enzymes
- Investigate effect of ubiquitination on substrate protein function



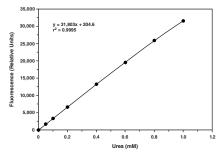
#### **Ubiquitinated Protein Capture Kit**

15979

- Demonstrate specific proteins are substrates for ubiquitin modification in vivo
- Capture and detect ubiquitinated proteins and free chains from cell lysates and tissue extracts by western blotting
- Identify and characterize ubiquitin-modified proteins by proteomic
- Investigate the role of ubiquitin in particular signaling pathways

#### Urea Fluorometric Assay Kit

- Measure urea in plasma, serum, and urine
- Assay 20 samples in duplicate
- Assay Range: 50 μM-1 mM
- Plate-based fluorometric measurement (ex 405-415 nm, em 470-480 nm)

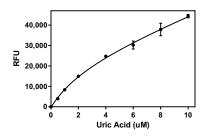


Uric Acid Assay Kit

700320 Vitellogenin (carp) ELISA Kit 10004993

10008883 WST-1 Cell Proliferation Assay Kit

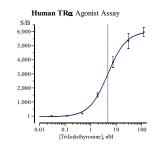
- Measure uric acid in serum, plasma, and urine
- Assay 40 samples in duplicate
- Assay Range: 0.5-10 μM
- Plate-based fluorometric measurement (ex 530-540 nm, em 585-595 nm)



#### VDR (human) Reporter Assay Kit

15742

- A nuclear receptor cell-based reporter assay
- See www.caymanchem.com for more details

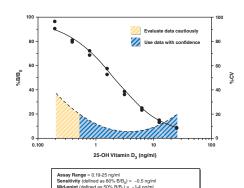




#### Vitamin D ELISA Kit

501050

- Measure vitamin D in plasma or serum
- Detect the metabolically stable forms 25(OH)D, and 25(OH)D,
- Assay 24 samples in triplicate or 36 samples in duplicate
- Measure vitamin D levels down to 0.5 ng/ml
- Incubation: 18 hours | Development: 60-90 minutes | Read: Colorimetric at 405-420 nm



ensitivity and mid-point were derived from the standa above. The standard was diluted with ELISA Buffer.

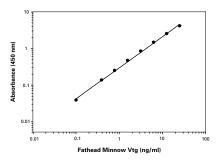
- An immunometric ELISA for analyzing vitellogenin (Vtg) in whole body homogenates and plasma samples
- See www.caymanchem.com for more details

#### bisense

#### Vitellogenin (fathead minnow) ELISA Kit

10006943

- An immunometric ELISA for analyzing vitellogenin (Vtg) in whole body homogenates
- See www.caymanchem.com for more details



bisense

#### Vitellogenin (medaka) ELISA Kit

10009223

- For the analysis of vitellogenin (Vtg) in whole body homogenates
- See www.caymanchem.com for more details

biôsense

#### Vitellogenin (rainbow trout) ELISA Kit

10004994

- An immunometric assay for the analysis of vitellogenin (Vtg)
- See www.caymanchem.com for more details

biosense

#### Vitellogenin (salmonid) Semi-Quantitative ELISA Kit

10009272

- A semi-quantitative assay for measuring vitellogenin (Vtg) from fish nlasma
- See www.caymanchem.com for more details

biosense

#### Vitellogenin (zebrafish) ELISA Kit

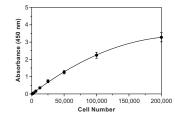
10004995

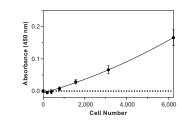
- An immunometric ELISA for analyzing vitellogenin (Vtg) in whole body homogenates and plasma samples
- See www.caymanchem.com for more details

biosense

#### • A colorimetric, non-radioactive assay for assessing cell viability and proliferation

- WST-1 is reduced to purple formazan by extracellular NADH from living cells
- Detect 1,000-200,000 cells per well
- Rapid assay; get results in 2-4 hours

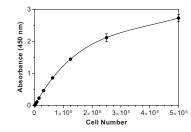




#### WST-8 Cell Proliferation Assay Kit

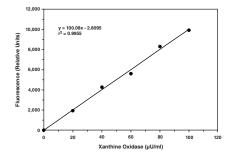
#### 10010199

- · A colorimetric, non-radioactive assay for assessing cell viability and proliferation
- WST-8 is reduced to purple formazan by extracellular NADH from living cells
- Detect 2,000-500,000 cells per well
- Rapid assay; get results in 2-4 hours

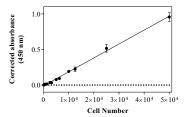


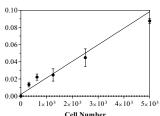
#### Xanthine Oxidase Fluorometric Assay Kit

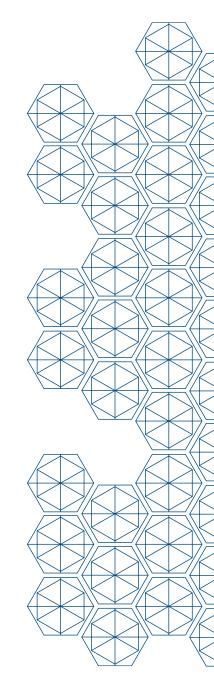
- Measure xanthine oxidase (XO) activity in plasma, serum, and tissue homogenates
- Assay 42 samples in duplicate
- Measure XO activity down to 10  $\mu$ U/ml
- Plate-based fluorometric measurement (ex 520-550 nm, em 585-595 nm)



- A colorimetric, non-radioactive assay for assessing cell viability and proliferation
- XTT is reduced to purple formazan by extracellular NADH from living cells
- Detect 500-100,000 cells per well
- Rapid assay; get results in about 4 hours







# **SAMPLE PREPARATION**

#### Angiotensin II Inhibitor Cocktail

9000681

- For storing blood samples containing angiotensin II
- Contains a mixture of inhibitors that prevent angiotensin II degradation





#### **CGRP Affinity Sorbent**

489009

- Immobilized mouse anti-CGRP antibody
- For use with the CGRP (rat) EIA Kit (Item No. 589001)



#### Cysteinyl Leukotriene Affinity Sorbent

400396

- Mouse monoclonal anti-LTC<sub>4</sub>/LTD<sub>4</sub>/LTE<sub>4</sub> antibody covalently bound to Sepharose 4B
- For affinity purification of cysteinyl LTs from complex matrices

#### **DNA Laddering Kit**

660990

- Maximal extraction of small DNA fragments from samples
- Minimal intact chromatin contamination

#### **Human Fibrinogen Affinity Sorbent**

401272

- Immobilized human fibrinogen
- For the removal of antibodies that bind unmodified fibrinogen
- For use with the Anti-Citrullinated Human Fibrinogen Assay Kit (mouse) (Item No. 501270)

#### Insulinase Inhibitor Cocktail

18429

- Allows insulin measurement even in hemolyzed samples
- · Contains insulinase inhibitors to protect sample integrity during collection



#### 8-Isoprostane Affinity Column

401111

- For purification of 8-isoprostane from urine and plasma
- Contains 0.5 ml of 8-isoprostane monoclonal antibody covalently bound to Sepharose beads
- Binding capacity is 5 ng of 8-isoprostane

#### 8-Isoprostane Affinity Purification Kit

501110

- Immunosorbent for the purification of 8-isoprostane from biological samples
- Employ a high affinity 8-isoprostane-specific monoclonal antibody

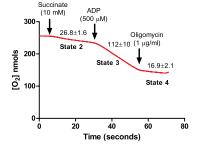
#### 8-Isoprostane Affinity Sorbent

• For purification of 8-isoprostane from urine and plasma

- Contains a 1 ml slurry of 8-isoprostane monoclonal antibody covalently bound to Sepharose beads in eicosanoid affinity column buffer
- Bind ~10 ng of 8-isoprostane/ml sorbent

#### MitoCheck® Mitochondrial (Tissue) Isolation Kit

- Isolate coupled mitochondria from fresh tissue samples
- Isolated mitochondria can be used with MitoCheck® Complex I-V Activity Assays Kits and Citrate Synthase Activity Assay Kit

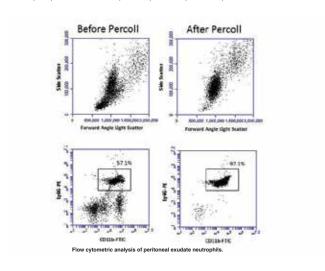


#### Neutrophil (mouse) Isolation Kit

601070

401113

- Isolate mouse neutrophils from peritoneal lavage or bone marrow
- Circumvents issues commonplace to isolating neutrophils from mouse blood
- Includes a Percoll density separation gradient
- Analyze purified neutrophils by flow cytometry



#### Nitrotyrosine Affinity Sorbent

389549

- Immobilized mouse anti-nitrotyrosine antibody
- For the capture of nitrated proteins from biological samples prior to SDS-PAGE

#### **Nuclear Extraction Kit**

10009277

- Isolate nuclear and cytoplasmic fractions from cell lysates and tissue homogenates
- Resulting isolates can be used in Transcription Factor Assay Kits
- Isolates can also be used for EMSA or western blotting
- Each kit provides reagents for 50 extractions from approximately 10 million cells

#### Sampling Tubes with BHT

• For storing oxidant-sensitive blood samples

• Contains the antioxidant butylated hydroxytoluene (BHT) (0.005% final concentration)



18534

10950

#### Sampling Tubes with BHT and Indomethacin

• For storing blood samples containing 8-isoprostanes and prostaglandin Contains the antioxidant butvlated hydroxytoluene (BHT) and the COX inhibitor indomethacin



#### Sampling Tubes with Indomethacin

10951

- For storing blood samples containing thromboxanes and prostaglandins
- Contains the COX inhibitor indomethacin (10 μM final concentration)





#### Sampling Tubes with PHMB

10949

- For storing cysteine protease-sensitive samples
- Contains the cysteine protease inhibitor polyhexamethylene biguanide (PHMB) (1 mM final concentration)





#### SPE Cartridges (C-18) (6 ml)

400020

- Solid phase extraction cartridges containing immobilized C-18
- For reverse-phase extraction of non-polar to moderately polar compounds
- Commonly used for the partial purification of eicosanoids

#### SPE Cartridges (Mixed Bed)

400021

- Solid phase extraction cartridges containing a mixture of non-polar and strong anion exchange chemistries
- For the partial purification of 11-dehydro TXB, from complex matrices

#### SPE Cartridges (Phenyl)

400048

- Solid phase extraction cartridges for aromatic and hydrophobic interactions
- For use with the Angiotensin II EIA Kit (Item No. 589301)



#### **SCREENING LIBRARIES**

#### Bio-active Lipid I Screening Library (96-Well)

- Consists of 11 plates with >820 compounds
- Includes prostaglandins, thromboxanes, cannabinoids, D-myo-inositol-phosphates, phosphatidylinositol-phosphates, sphingolipids, ceramide derivatives, inhibitors, and receptor agonists and antagonists

#### Bio-active Lipid II Screening Library (96-Well)

- Consists of 3 plates with >190 compounds
- Includes prostaglandins, isoprostanes, thromboxanes, leukotrienes, lipoxins, and several other complex lipid mediators

#### **Epigenetics Screening Library (96-Well)**

- Contains >140 small molecules
- Includes compounds that modulate the activity of methyltransferases, demethylases, HATs, HDACs, and acetylated lysine reader proteins

#### Fatty Acid Screening Library (96-Well)

- Contains > 75 compounds
- Includes arachidonic, linoleic, eicosapentaenoic, stearic, α- and γ-linolenic, oleic, and palmitic acids

#### Kinase Screening Library (96-Well)

10505

10506

10507

11076

10504

- Consists of 2 plates with >140 specific and non-specific kinase inhibitors
- Includes dual specific threonine/tyrosine kinase inhibitors, selective MAPK inhibitors, p38 MAPK inhibitors, and PI3K inhibitors

#### Prostaglandin Screening Library I (96-Well) 10501

- Contains > 70 compounds
- Includes the "F-series" prostaglandins and isoprostanes

#### 10502 Prostaglandin Screening Library II (96-Well)

- Contains >60 compounds
- Includes the "D and E-series" prostaglandins

#### Prostaglandin Screening Library III (96-Well) 10503

- Contains >75 compounds
- Includes the "A and J-series" prostaglandins

#### Stem Cell Small Molecule Screening Library (96-Well)

9001827

- Contains >140 small molecules in two 96-well plates
- Includes compounds that induce differentiation, maintain self-renewal and proliferation, or improve the reprogramming efficiency of various stem cell populations

\*The composition of the libraries will always vary somewhat depending upon our inventory.



# ALPHABETICAL INDEX

$\triangle$	
7-AAD/CFSE Cell-Mediated Cytotoxicity Assay Kit	15
Acid Phosphatase Colorimetric Assay Kit	
Aconitase Assay Kit	
AcSDKP EIA Kit	
Acylated Ghrelin (dog) Easy Sampling EIA Kit	
Acylated Ghrelin (human) Easy Sampling EIA Kit	
Acylated Ghrelin (human) Express EIA Kit	
Acylated Ghrelin (mouse, rat) Easy Sampling EIA Kit	
Acylated Ghrelin (mouse, rat) Express EIA Kit	
Acylated Ghrelin (pig) EIA Kit	
Adipogenesis Assay Kit	
Adipolysis Assay Kit	
Adiponectin (human) EIA Kit	
Adiponectin (human) EIA Kit (HS)	
Adiponectin (mouse) EIA Kit	
AhR (human) Assay Kit	
Alanine Transaminase Colorimetric Activity Assay Kit	
Aldehyde Dehydrogenase Activity Assay Kit	
Aldosterone ELISA Kit	
Angiotensin II EIA Kit	
Angiotensin II EIA Nit	
Angiotensin'il lillibitor Cocktail	
Anti-Citrullinated Human Fibrinogen Assay Kit (mouse)	
Anti-Ovalbumin IgE (mouse) ELISA Kit Anti-Ovalbumin IgG, (mouse) ELISA Kit	
Antioxidant Assay KitApoA1 (human) ELISA Kit	
AR (human) Reporter Assay Kit	
Arginine Vasopressin ELISA Kit	
Ascorbate Assay Kit	
ASP ELISA Kit	
Atriopeptin (rat) EIA Kit	
Autophagy/Cytotoxicity Dual Staining Kit	
Autophlux Kit	
Autotaxin Inhibitor Screening Assay Kit	
Autotaxiii iiiiibitoi Scieeliilig Assay Nt	17
В	
BAZ2B bromodomain TR-FRET Assay Kit	19
Bio-active Lipid I Screening Library (96-Well)	
Bio-active Lipid II Screening Library (96-Well)	
BRD2 bromodomain 1 TR-FRET Assay Kit	
BRD2 bromodomain 2 TR-FRET Assay Kit	
BRD2 bromodomains 1 and 2 TR-FRET Assay Kit	
BRD3 bromodomain 1 TR-FRET Assay Kit	
BRD3 bromodomain 2 TR-FRET Assay Kit	
BRD3 bromodomains 1 and 2 TR-FRET Assay Kit	
BRD4 bromodomain 1 TR-FRET Assay Kit	
BRD4 bromodomain 2 TR-FRET Assay Kit	
BRD4 bromodomains 1 and 2 TR-FRET Assay Kit	
BRDT bromodomain 1 TR-FRET Assay Kit	
BRG1 bromodomain TR-FRET Assay Kit	
BRM bromodomain TR-FRET Assay Kit	
,	
Calcium Assay Kit	21
CAR1 (human) Reporter Assay Kit	
CAR3 (human) Reporter Assay Kit	
Caspase-3 Fluorescence Assay Kit	
Catalase Assay Kit	
Catalase Assay Kit (without Hydrogen Peroxide)	22
Cayman Practice ELISA Kit	22
CBP bromodomain TR-FRET Assay Kit	

Cell Cycle Phase Determination Kit	.23
CFSE Cell Division Assay Kit	.23
CGRP Affinity Sorbent	
CGRP (human) EIA Kit	
CGRP (rat) EIA Kit	
Cholesterol Cell-Based Detection Assay Kit	
Cholesterol Fluorometric Assay Kit	
Cholesterol Uptake Cell-Based Assay KitAritimusessessessessessessessessessessessesses	
Corticosterone ELISA Kit	
Cortisol ELISA Kit	
Cortisol Express ELISA Kit	
COX Activity Assay Kit	
COX Colorimetric Inhibitor Screening Assay Kit	
COX (ovine) Colorimetric Inhibitor Screening Assay Kit	
COX Fluorescent Activity Assay Kit	
COX Fluorescent Inhibitor Screening Assay Kit	
COX (ovine) Inhibitor Screening Assay Kit	
COX (ovine/human) Inhibitor Screening Assay Kit	
COX-1 (human) Inhibitor Screening Assay Kit	
COX-1 (ovine) Inhibitor Screening Assay KitCOX-2 (human) Inhibitor Screening Assay Kit	
COX-2 (numar) Inhibitor Screening Assay Kit	
cPLA, Assay Kit	
C-Reactive Protein (human) ELISA Kit	
Creatinine (serum) Colorimetric Assay Kit	
Creatinine (urinary) Colorimetric Assay Kit	
CREB (Phospho-Ser <sup>133</sup> ) Transcription Factor Assay Kit	
Cyclic AMP ELISA Kit	
Cyclic AMP ELISA Kit (without Acetic Anhydride)	
Cyclic AMP Select ELISA Kit	
Cyclic GMP ELISA Kit	
Cyclic GMP ELISA Kit (without Acetic Anhydride)	
CYP1A1/2 Induction Reporter Assay KitCYP2B6 Induction Reporter Assay Kit	
CYP2C9 Induction Reporter Assay KitCYP2C9 Induction Reporter Assay Kit	
Cysteinyl Leukotriene Affinity Sorbent	
Cysteinyl Leukotriene ELISA Kit	
Cysteinyl Leukotriene Express ELISA Kit	
11-dehydro Thromboxane B <sub>2</sub> ELISA Kit - Monoclonal	73
DHN-MA EIA Kit	
13,14-dihydro-15-keto Prostaglandin F <sub>2a</sub> ELISA Kit	.64
DNA Laddering Kit	
DNA Methylation ELISA Kit	
DNA/RNA Oxidative Damage ELISA Kit	.29
DPP (IV) Inhibitor Screening Assay Kit	
DUB Activity Assay Kit	.30
<del>_</del>	
_	
 Endothelin ELISA Kit	.30
Enterolactone ELISA Kit	.30
EP <sub>2</sub> Receptor (rat) Reporter Assay Kit	.30
EP4 Receptor (rat) Activation Assay Kit (cAMP)	.30
EP <sub>4</sub> Receptor (rat) Reporter Assay Kit	.31
Epigenetics Screening Library (96-Well)	
EpiTitan <sup>TM</sup> Histone Peptide Array	
ER (human) Reporter Assays Panel	
ERα (human) Reporter Assay Kit	
ERβ (human) Reporter Assay Kit	
ERRα (human) Reporter Assay Kit ERRγ (human) Reporter Assay Kit	
Estradiol ELISA Kit	
Estrial FLICA I/it	22

F	
FAAH Inhibitor Screening Assay Kit	32
FABP4 (human) EIA Kit	
FABP4 Inhibitor/Ligand Screening Assay Kit	
Fatty Acid Screening Library (96-Well)	
Fluprostenol ELISA Kit	
Follicle-Stimulating Hormone ELISA Kit	33
Free Fatty Acid Fluorometric Assay Kit	33
FXR (human) Reporter Assay Kit	33
G	
GFAP (human) EIA Kit	33
Acylated Ghrelin (dog) Easy Sampling EIA Kit	
Acylated Ghrelin (human) Easy Sampling EIA Kit	
Acylated Ghrelin (human) Express EIA Kit	
Acylated Ghrelin (mouse, rat) Easy Sampling EIA K	
Acylated Ghrelin (mouse, rat) Express EIA Kit	
Acylated Ghrelin (pig) EIA Kit	
Unacylated Ghrelin (dog) Easy Sampling EIA Kit	34
Unacylated Ghrelin (human) Easy Sampling EIA Ki	t35
Unacylated Ghrelin (mouse, rat) Easy Sampling El	
Unacylated Ghrelin (pig) EIA Kit	
Unacylated Ghrelin (human) Express EIA Kit	
Unacylated Ghrelin (mouse, rat) Express EIA Kit	
GLP SAM-Screener™ Assay Kit	
Glucose Colorimetric Assay Kit	
Glucose-6-Phosphate Dehydrogenase Activity Ass	'
Glucose-6-Phosphate Fluorometric Assay Kit Glucose Uptake Cell-Based Assay Kit	
Glutaredoxin Fluorescent Activity Assay Kit	
Glutathione Assay Kit	
Glutathione Cell-Based Detection Kit (Blue Fluore	
Glutathione Peroxidase Assay Kit	,
Glutathione Reductase Assay Kit	
Glutathione S-Transferase Assay Kit	
S-Glutathionylated Protein Detection Kit	37
Glycerol Cell-Based Assay Kit	37
Glycerol Colorimetric Assay Kit	
Glycogen Assay Kit	
Glycolysis Cell-Based Assay Kit	
Glyco-SPOT DNA Repair Assay Kit	
Goat Anti-Human Fluorophore Selection Kit	
Goat Anti-Robbit Fluorophore Selection Kit	
Goat Anti-Rabbit Fluorophore Selection Kit GR (human) Reporter Assay Kit	30 37
Growth Hormone (rat) EIA Kit	
Growth From the traty EIA Nit	
1.1	
HAT Inhibitor Screening Assay Kit	
HDAC Cell-Based Activity Assay Kit	
HDAC Fluorometric Activity Assay Kit	
HDAC1 Inhibitor Screening Assay Kit	
HDAC8 Inhibitor Screening Assay Kit	
Hemoglobin Colorimetric Assay Kit	
15(S)-HETE ELISA Kit HIF-1α Transcription Factor Assay Kit	
His-Tag Detection ELISA Kit	
Histamine EIA Kit	
Human Fibrinogen Affinity Sorbent	
Human Serum Albumin ELISA Kit	
Human Therapeutic IgG1 ELISA Kit	
Human Therapeutic IgG2 ELISA Kit	
Human Therapeutic IgG4 ELISA Kit	
Hydrogen Peroxide Cell-Based Assay Kit	
β-Hydroxybutyrate (Ketone Body) Colorimetric As	ssay Kit42
β-Hydroxybutyrate (Ketone Body) Fluorometric A	ssay Kit42
20-Hydroxyecdysone EIA Kit	
8-Hydroxyguanosine Monoclonal FLISA Kit	42

nsulin (rat) EIA Kit	42
nsulinase Inhibitor Cocktail	77
nterleukin-1α (human) ELISA Kit	
nterleukin-1β (human) ELISA Kit	
nterleukin-6 (human) ELISA Kit	
nterleukin-6 (mouse) ELISA Kit	
3-Isoprostane Affinity Column 3-Isoprostane Affinity Purification Kit	
3-Isoprostane Affinity Sorbent	
B-Isoprostane ELISA Kit	
3-Isoprostane Express ELISA Kit	
1	
J	
IC-1 Mitochondrial Membrane Potential Assay Kit	
IMJD2A Tudor Domains TR-FRET Assay Kit	44
6-keto Prostaglandin F <sub>1α</sub> ELISA Kit	
11-keto Testosterone ELISA Kit	
Kinase Screening Library (96-Well)	/8
 _3MBTL1 MBT Domains TR-FRET Assay Kit	44
D-Lactate Assay Kit	
-Lactate Assay Kit	
_atanoprost ELISA Kit	
_C3 Interact Kit	45
_DH Cytotoxicity Assay Kit	
_DL Uptake Cell-Based Assay Kit	
Leptin (human) EIA Kit	
eptin (mouse/rat) EIA Kit	
eptin Receptor (human) EIA Kit	
Leukotriene $\rm B_4$ ELISA Kiteukotriene $\rm B_4$ Express ELISA Kit	4040 4.
Leukotriene C <sub>4</sub> ELISA Kit	
Leukotriene E <sub>4</sub> ELISA Kit	
ipase Activity Assay Kit	
Lipid Droplets Fluorescence Assay Kit	
_ipid Hydroperoxide (LPO) Assay Kit	
Lipid Hydroperoxide (LPO) Assay Kit (96 well)	
Lipoxygenase Inhibitor Screening Assay Kit	
Live Cell Multiplex Assay Kit	
LSD1 Inhibitor Screening Assay Kituteinizing Hormone ELISA Kit	
XR (human) Reporter Assays Panel	
XRα (human) Reporter Assay Kit	
XRβ (human) Reporter Assay Kit	
× 1	
$\vee$	
Malachite Green Phosphate Assay Kit	
Melanocortin-3 Receptor Reporter Assay Kit	
Melanocortin-4 Receptor Reporter Assay Kit	
2-Methoxyestradiol ELISA Kit	
Methyltransferase Colorimetric Assay Kit	
Methyltransferase Fluorometric Assay KitMitoCheck® Citrate Synthase Activity Assay Kit	
MitoCheck® Complex   Activity Assay Kit	
MitoCheck® Complex I Activity Assay Kit	
MitoCheck® Complex II/III Activity Assay Kit	
MitoCheck® Complex IV Activity Assay Kit	
MitoCheck® Complex V Activity Assay Kit	50
MitoCheck® Mitochondrial (Tissue) Isolation Kit	
MitoCheck® Mitochondrial OCR Assay Kit	
MLL1 SAM-Screener <sup>™</sup> Assay Kit	50

Mouse Anti-Type II Collagen IgG Assay Kit (bovine)	
Mouse IgG ELISA Kit	
MR (human) Reporter Assay Kit	
MTT Cell Proliferation Assay Kit	
Multi-Drug Resistance Assay Kit (Calcein AM)	
Myeloperoxidase Chlorination Fluorometric Assay Kit	
Myeloperoxidase Inhibitor Screening Assay Kit	
Myeloperoxidase Peroxidation Fluorometric Assay Kit	
NAD/NADH Cell-Based Assay Kit	52
Neutrophil Elastase Activity Assay Kit	
Neutrophil Extracellular Trap (MPO-Histone) ELISA Kit	
Neutrophil Extracellular Trap (NET) Assay Kit	
Neutrophil (mouse) Isolation Kit	
Neutrophil/Monocyte Respiratory Burst Assay Kit	
Neutrophil Myeloperoxidase Activity Assay Kit	
NF-кВ (human p50) Transcription Factor Assay KitNF-кВ (p65) Transcription Factor Assay Kit	
Nitrate/Nitrite Colorimetric Assay Kit	
Nitrate/Nitrite Colorimetric Assay Kit (LDH method)	
Nitrate/Nitrite Fluorometric Assay Kit	
S-Nitrosylated Protein Detection Kit	.54
Nitrotyrosine Affinity Sorbent	
NOS Activity Assay Kit	
Nrf2 (human) Reporter Assay Kit	
Nrf2 Transcription Factor Assay Kit	
Nucleal Extraction Nit	. / /
$\bigcup$	
Oct4 Transcription Factor Assay Kit	
Orexin 1 Receptor Reporter Assay Kit Orexin 2 Receptor Reporter Assay Kit	
Anti-Ovalbumin IgE (mouse) ELISA Kit	
Anti-Ovalbumin IgG, (mouse) ELISA Kit	
Oxygen Consumption Rate Assay Kit (MitoXpress®-Xtra HS Method)	
Oxygen Consumption/Glycolysis Dual Assay Kit	
Oxygen Consumption/MitoMembrane Potential Dual Assay Kit	
Oxytocin ELISA Kit	.56
P	
p53 Transcription Factor Assay Kit	.56
PAD4 Autoantibody ELISA Kit	.56
PAD4 Inhibitor Screening Assay Kit (Ammonia)	
PAF Acetylhydrolase Assay Kit	.56
PAF Acetylhydrolase Inhibitor Screening Assay Kit	
tetranor-PGDM ELISA KitPGR (human) Reporter Assay Kit	
Phagocytosis Assay Kit (IgG FITC)	
Phagocytosis Assay Kit (IgG PE)	
17-phenyl trinor Prostaglandin F <sub>2α</sub> ELISA Kit	.64
Phosphatidylcholine Colorimetric Assay Kit	.57
cPLA <sub>2</sub> Assay Kit	
sPLA <sub>2</sub> Assay Kit	
sPLA <sub>2</sub> (human Type IIA) ELISA Kit	.58
sPLA <sub>2</sub> (Type V) Inhibitor Screening Assay Kit	.JO 58
PPARα (human) Reporter Assay Kit	
PPARα (mouse) Reporter Assay Kit	
PPARα (mouse) Reporter Assay Kit	.58
PPARα (rat) Reporter Assay Kit PPARα Transcription Factor Assay Kit	.58 .58 .59
PPARα (rat) Reporter Assay Kit PPARα Transcription Factor Assay Kit PPARα, δ, γ Complete Transcription Factor Assay Kit	.58 .58 .59
PPARα (rat) Reporter Assay Kit	.58 .58 .59 .59
PPARα (rat) Reporter Assay Kit	.58 .58 .59 .59 .59
PPARα (rat) Reporter Assay Kit	.58 .59 .59 .59 .59
PPARα (rat) Reporter Assay Kit	.58 .59 .59 .59 .59 .59

PPARγ (mouse/rat) Reporter Assay Kit	60
PPARy Transcription Factor Assay Kit	
Pregnane X (human) Receptor Assay Kit	
Pregnane X (rat) Receptor Assay Kit	
Progesterone ELISA Kit	
Prolactin (human) ELISA Kit	
Prolactin (rat) EIA Kit	
Prostaglandin D Synthase (hematopoietic-type) FP-Based	OI
	/ 1
Inhibitor Screening Assay Kit - Green	/
Prostaglandin D <sub>2</sub> ELISA Kit	61
Prostaglandin D <sub>2</sub> Express ELISA Kit	61
Prostaglandin D <sub>2</sub> -MOX ELISA Kit	
Prostaglandin D <sub>2</sub> -MOX Express ELISA Kit	62
Prostaglandin E Metabolite ELISA Kit	
Prostaglandin E <sub>2</sub> ELISA Kit - Monoclonal	
Prostaglandin E <sub>2</sub> Express ELISA Kit	
6-keto Prostaglandin F <sub>1a</sub> ELISA Kit	63
Prostaglandin F., ELISA Kit	63
11β-Prostaglandin $F_{2\alpha}$ ELISA Kit	63
13.14-dihvdro-15-keto Prostaglandin F. ELISA Kit	64
17-phenyl trinor Prostaglandin E FLISA Kit	64
17-phenyl trinor Prostaglandin F $_{2\alpha}$ ELISA Kit Prostaglandin I Metabolite ELISA Kit	64
Prostaglandin Screening ELISA Kit	6/1
Prostaglandin Screening Library I (96-Well)	
Prostaglandin Screening Library II (96-Well)	
Prostaglandin Screening Library III (96-Well)	
20S Proteasome Assay Kit	
Protein Carbonyl Colorimetric Assay Kit	
Protein Determination Kit	
Protein Synthesis Assay Kit	
Pyruvate Assay Kit	65
R	
Rabbit Anti-Goat Fluorophore Selection Kit	65
RAR (human) Reporter Assays Panel	65
RARα (human) Reporter Assay Kit	65
RARβ (human) Reporter Assay Kit	66
RARγ (human) Reporter Assay Kit	
C-Reactive Protein (human) ELISA Kit	
Renin Inhibitor Screening Assay Kit	
Resistin (human) EIA Kit	
Resistin (rat) EIA Kit	
Resolvin D1 ELISA Kit	
Resolvin D2 ELISA Kit	
RORγ (human) Reporter Assay Kit	
RXR (human) Reporter Assays Panel	67
RXRα (human) Reporter Assay Kit	67
RXRβ (human) Reporter Assay Kit	67
RXRy (human) Reporter Assay Kit	67
C	
$\supset$	
Sampling Tubes with BHT	78
Sampling Tubes with BHT and Indomethacin	
Sampling Tubes with Indomethacin	
Sampling Tubes with PHMB	
Secreted Alkaline Phosphatase Reporter Gene	, 0
Assay Kit (Luminescence)	68
Semi-Quantitative Biomarker ELISA Component Kit (anti-mouse)	
Semi-Quantitative Biomarker ELISA Component Kit (anti-rabbit)	
SET7/9 SAM-Screener™ Assay Kit	
SIRT1 Direct Fluorescent Screening Assay Kit	
SIRT1 FRET-Based Screening Assay Kit	
SIRT2 Direct Fluorescent Screening Assay Kit	
SIRT3 Direct Fluorescent Screening Assay Kit	
Soluble Epoxide Hydrolase Cell-Based Assay Kit	69
Soluble Epoxide Hydrolase Inhibitor Screening Assay Kit	
SPE Cartridges (C-18) (6 ml)	
SPE Cartridges (Mixed Bed)	
01 2 941 41 46 600 (1 11) 404 204 (1 11)	

SPE Cartridges (Phenyl)	
Sphingomyelin Colorimetric Assay Kit	
sPLA <sub>2</sub> Assay Kit	
sPLA <sub>2</sub> (human Type IIA) ELISA Kit	
sPLA <sub>2</sub> (Type V) Inhibitor Screening Assay Kit	
SREBP-1 Transcription Factor Assay Kit SREBP-2 Cell-Based Translocation Assay Kit	
SREBP-2 Transcription Factor Assay Kit	
STAT-8-Isoprostane ELISA Kit	
Steatosis Colorimetric Assay Kit	
Stem Cell Small Molecule Screening Library (96-W	
Substance P ELISA Kit	
Sulfenylated Protein Cell-Based Detection Kit	
Superoxide Dismutase Assay Kit	
SureLight® APC Conjugation Kit 1 mg lgG	7
SureLight® APC Conjugation Kit 2 mg lgG	
SureLight® R-PE Conjugation Kit	7
T	
TAF1 bromodomain 1 TR-FRET Assay Kit	7
TAF1 bromodomains 1 and 2 TR-FRET Assay Kit	
TBARS Assay Kit	
TBARS (TCA Method) Assay Kit	
Testosterone ELISA Kit	
11-keto Testosterone ELISA Kit	
Thiol Detection Assay Kit	
Thioredoxin Activity Fluorescent V2 Assay Kit	
Thioredoxin Reductase Colorimetric Assay Kit	
Thioredoxin Reductase Fluorescent Activity Assay	Kit7
Thioredoxin/Thioredoxin Reductase Mammalian A	
Thromboxane B <sub>2</sub> ELISA Kit	
Thromboxane B <sub>2</sub> Express ELISA Kit - Monoclonal	
11-dehydro Thromboxane B <sub>2</sub> ELISA Kit - Monoclo	
Thyroid-Stimulating Hormone ELISA KitTMRE Mitochondrial Membrane Potential Assay K	
TNF-α (human) ELISA Kit	
TNF-a (mouse) ELISA Kit	
TR (human) Reporter Assays Panel	
TRα (human) Reporter Assay Kit	
TRβ (human) Reporter Assay Kit	
Triglyceride Colorimetric Assay Kit	
	_
Ubiquitin Interact Kit	
Ubiquitin Link KitUbiquitinated Protein Capture Kit	
Unacylated Ghrelin (dog) Easy Sampling EIA Kit	
Unacylated Ghrelin (dog) Easy Sampling EIA Rit  Unacylated Ghrelin (human) Easy Sampling EIA Rit	
Unacylated Ghrelin (mouse, rat) Easy Sampling EIA	
Unacylated Ghrelin (pig) EIA Kit	
Unacylated Ghrelin (human) Express EIA Kit	
Unacylated Ghrelin (mouse, rat) Express EIA Kit	
Urea Fluorometric Assay Kit	7
Uric Acid Assay Kit	7
V VDD (human) Danartar Assaul/it	-
VDR (human) Reporter Assay KitVitamin D ELISA Kit	
Vitellogenin (carp) ELISA Kit	
Vitellogenin (fathead minnow) ELISA Kit	
Vitellogenin (medaka) ELISA KitVitellogenin (medaka) ELISA Kit	
Vitellogenin (rainbow trout) ELISA Kit	
Vitellogenin (salmonid) Semi-Quantitative ELISA K	
Vitellogenin (zebrafish) ELISA Kit	

$\bigvee$	
WST-1 Cell Proliferation Assay Kit	
X	
Xanthine Oxidase Fluorometric Assay Kit	





# ORDERS & TECHNICAL SUPPORT

USA & Canada www.caymanchem.com Toll Free (800) 364-9897

International Distributors www.caymanchem.com/distributors

Technical Support www.caymanchem.com/tech Toll Free (888) 526-5351

