

Research Tools for Lipid Analysis by Mass Spectrometry



Mass spectrometry is a powerful technique to characterize the lipid profile in biological samples. High-quality lipid reference standards are essential during mass spectrometry-based lipid analysis to produce accurate, reliable, and reproducible results. Cayman has been a trusted source of lipid research standards for over 45 years. Backed by decades of experience in lipid synthesis, purification, and characterization, our high-purity lipid standards have helped researchers explore fundamental mechanisms of biology, identify biomarkers, and develop new therapies.

Explore our Lipid Standards & Custom Research Services

- High-purity, qualitative research standards
- Stable isotope-labeled standards
- Ready-to-use MaxSpec[®] analytical standard mixtures
- Untargeted lipidomics & targeted lipid analysis
- Custom lipid synthesis



New

Deuterated Lipidomics MaxSpec[®] Mixture

Item No. 40974

- Contains 15 lipid class-specific internal standards for signal normalization and semiquantification
- Enables absolute quantification when paired with authentic unlabeled standards
- Confirms MS/MS and retention time method performance during method development and validation

New

MaxSpec[®] Reference Plasma (Human)

Item No. 42573

A reference material for the analysis of lipids or metabolites

Use for:

- Optimizing instrument parameters
- Monitoring instrument precision
- Evaluating recovery from extraction
- Validating overall method accuracy

MaxSpec[®] Mixtures for LC-MS

Ready-to-Use Quantitative Standards

Our MaxSpec[®] mixtures streamline workflows by reducing the effort and variability of mixing individual internal and calibration standards.

MaxSpec[®] Quality Guarantee

- ✓ Headspace purged with argon
- ✓ At least 1 ml of solution per ampule
- ✓ Packaged in glass ampoules
- ✓ Custom mixtures also available
- ✓ Designed for direct snap-and-inject use

Fatty Acyls

Prostaglandins

Unlabeled Standards

Item No.	Product Name
14010	Prostaglandin E ₂
16010	Prostaglandin F _{2α}
18500	Prostaglandin J ₂
12010	Prostaglandin D ₂

Popular matched pair items shown, over 435 prostaglandins, intermediates, and metabolites available online

Labeled Standards

Item No.	Product Name
314010	Prostaglandin E ₂ -d ₄
316010	Prostaglandin F _{2α} -d ₄
19345	Prostaglandin J ₂ -d ₄
312010	Prostaglandin D ₂ -d ₄

MaxSpec® Mixtures for Prostaglandins

Item No.	Product Name
19300	Primary Prostaglandin Metabolite MaxSpec® LC-MS Mixture*

Item No.	Product Name
19422	Secondary Prostaglandin Metabolite MaxSpec® LC-MS Mixture

* Also Available: Deuterated Primary Prostaglandin Metabolite MaxSpec® LC-MS Mixture (Item No. 19304)

Hydroxy/Hydroperoxy/Epoxy/Oxo Eicosanoids

Unlabeled Standards

Item No.	Product Name
34230	5(S)-HETE
34570	12(S)-HETE
34720	15(S)-HETE
90030	20-HETE
50511	(±)11(12)-EET
50651	(±)14(15)-EET
51651	(±)14(15)-DiHET
34250	5-OxoETE

Popular matched pair items shown, over 170 hydroxy/hydroperoxy/epoxy/oxo eicosanoids available online

Labeled Standards

Item No.	Product Name
334230	5(S)-HETE-d ₈
334570	12(S)-HETE-d ₈
334720	15(S)-HETE-d ₈
390030	20-HETE-d ₆
10006413	(±)11(12)-EET-d ₁₁
10006410	(±)14(15)-EET-d ₁₁
10008040	(±)14(15)-DiHET-d ₁₁
334250	5-OxoETE-d ₇

Thromboxanes & Leukotrienes

Unlabeled Standards

Item No.	Product Name
19030	Thromboxane B ₂
19500	11-dehydro Thromboxane B ₂
20110	Leukotriene B ₄
20210	Leukotriene C ₄

Popular matched pair items shown, over 20 thromboxanes and 40 leukotrienes available online

Labeled Standards

Item No.	Product Name
319030	Thromboxane B ₂ -d ₄
319500	11-dehydro Thromboxane B ₂ -d ₄
320110	Leukotriene B ₄ -d ₄
10006198	Leukotriene C ₄ -d ₅

MaxSpec® Mixtures for Arachidonic Acid Metabolites

Item No.	Product Name
19101	Primary COX and LOX MaxSpec® LC-MS Mixture*

Item No.	Product Name
20665	Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture

Item No.	Product Name
20666	Arachidonic Acid Oxylipin MaxSpec® LC-MS Mixture*

* Also available: Deuterated Primary COX and LOX MaxSpec® LC-MS Mixture (Item No. 19228) and Deuterated Arachidonic Acid Oxylipin MaxSpec® LC-MS Mixture (Item No. 20668)

Octadecanoids

Unlabeled Standards

Item No.	Product Name
38610	13(S)-HODE
38410	9(S)-HODE
52400	(±)9(10)-EpOME
53400	(±)9(10)-DiHOME
52450	(±)12(13)-EpOME
10009832	(±)12(13)-DiHOME

Popular matched pair items shown, over 250 octadecanoids available online

Labeled Standards

Item No.	Product Name
338610	13(S)-HODE-d ₄
338410	9(S)-HODE-d ₄
10009995	(±)9(10)-EpOME-d ₄
10009993	(±)9(10)-DiHOME-d ₄
10009996	(±)12(13)-EpOME-d ₄
10009994	(±)12(13)-DiHOME-d ₄

MaxSpec® Mixtures for Octadecanoids

Item No.	Product Name
20794	Linoleic Acid Oxylipins MaxSpec® LC-MS Mixture*

Item No.	Product Name
22638	ALA and GLA Oxylipin MaxSpec® LC-MS Mixture

Item No.	Product Name
22281	Dihomo-γ-Linolenic Acid Metabolite MaxSpec® LC-MS Mixture

* Also available: Deuterated Linoleic Acid Oxylipins MaxSpec® LC-MS Mixture (Item No. 20795)

Specialized Pro-Resolving Mediators

Docosanoids

Unlabeled Standards

Item No.	Product Name
13060	17(R)-Resolvin D1
10009799	17(S)-HDHA
10878	Maresin 1
10010390	Protectin D1
10012554	Resolvin D1
13834	Resolvin D3

Popular matched pair items shown, over 145 docosanoids available online

Labeled Standards

Item No.	Product Name
11183	17(R)-Resolvin D1-d ₅
29797	17(S)-HDHA-d ₅
21823	Maresin 1-d ₅
21625	Protectin D1-d ₅
11182	Resolvin D1-d ₅
19512	Resolvin D3-d ₅

Lipoxins & E-Series Resolvins

Unlabeled Standards

Item No.	Product Name
90410	Lipoxin A ₄
10007848	Resolvin E1

Popular matched pair items shown, over 10 lipoxins and resolvin Es available online

Labeled Standards

Item No.	Product Name
10007737	Lipoxin A ₄ -d ₅
10009854	Resolvin E1-d ₄

MaxSpec® Mixtures for Specialized Pro-Resolving Mediators

Item No.	Product Name
19412	Lipoxin MaxSpec® LC-MS Mixture

Item No.	Product Name
18702	SPM D-series MaxSpec® LC-MS Mixture

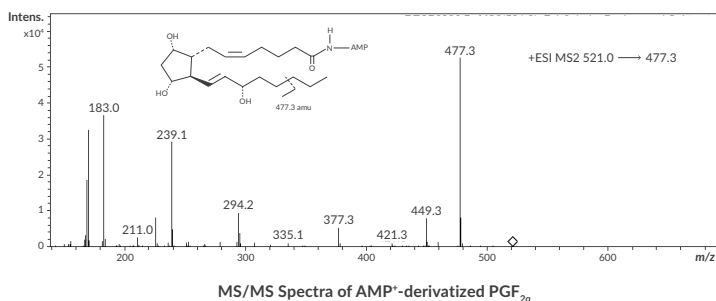
Item No.	Product Name
19417	SPM E-series MaxSpec® LC-MS Mixture

AMP⁺ MaxSpec[®] Kit

Item No. 710000

A novel method to increase MS sensitivity for detection of lipids

- AMP⁺ can be coupled quantitatively to free carboxylic acids through an amide bond linkage
- Derivatized product shows improved electrospray ionization and improved MS sensitivity
- Stable to MS/MS fragmentation; isobaric derivatives yield unique fragment ions



Fatty Acids

Unlabeled Standards

Item No.	Product Name
13351	Myristic Acid; C14:0
10006627	Palmitic Acid; C16:0
10011298	Stearic Acid; C18:0
90260	Oleic Acid; C18:1
90150	Linoleic Acid; C18:2
90210	α-Linolenic Acid; C18:3
90230	Dihomo-γ-Linolenic Acid; C20:3
90010	Arachidonic Acid; C20:4
90110	Eicosapentaenoic Acid; C20:5
90310	Docosahexaenoic Acid; C22:6
13353	Lignoceric Acid; C24:0

Popular matched pair items shown, over 340 fatty acids available online

Labeled Standards

Item No.	Product Name
9003317	Myristic Acid-d ₂₇ ; C14:0-d ₂₇
16497	Palmitic Acid-d ₃₁ ; C16:0-d ₃₁
9000455	Stearic Acid-d ₅ ; C18:0-d ₅
9000432	Oleic Acid-d ₁₇ ; C18:1-d ₁₇
9002193	Linoleic Acid-d ₁₁ ; C18:2-d ₁₁
9002286	α-Linolenic Acid-d ₅ ; C18:3-d ₅
10458	Dihomo-γ-Linolenic Acid-d ₆ ; C20:3-d ₆
390010	Arachidonic Acid-d ₈ ; C20:4-d ₈
10005056	Eicosapentaenoic Acid-d ₅ ; C20:5-d ₅
10005057	Docosahexaenoic Acid-d ₅ ; C22:6-d ₅
9003320	Lignoceric Acid-d ₄₇ ; C24:0-d ₄₇

MaxSpec[®] Mixtures for Fatty Acids

Item No.	Product Name
17941	Polyunsaturated Fatty Acid MaxSpec [®] LC-MS Mixture
17942	Saturated/Monounsaturated Fatty Acid MaxSpec [®] LC-MS Mixture

Item No.	Product Name
18824	Short-Chain Fatty Acid MaxSpec [®] LC-MS Mixture
19227	Odd-Chain Fatty Acid MaxSpec [®] LC-MS Mixture

Edible Oil Mixtures

Item No.	Product Name
29365	AOCS Canola Oil Fatty Acid Methyl Ester Standard Mixture
28665	AOCS Long-Chain Fatty Acid Methyl Ester Standard Mixture 1
28666	AOCS Long-Chain Fatty Acid Methyl Ester Standard Mixture 2
29366	AOCS Long-Chain Fatty Acid Methyl Ester Standard Mixture 3
28668	AOCS Long-Chain Fatty Acid Methyl Ester Standard Mixture 4
28669	AOCS Medium- and Long-Chain Fatty Acid Methyl Ester Standard Mixture 1
28670	AOCS Medium- and Long-Chain Fatty Acid Methyl Ester Standard Mixture 2

Lipidomics & Lipid Analysis Services

Cayman's lipid analysis laboratories are equipped with state-of-the-art chromatography/mass spectrometry systems and data analysis software, enabling us to offer lipidomics and lipid analysis services to meet your specific project needs.

Learn more at www.caymanchem.com/lipidomics

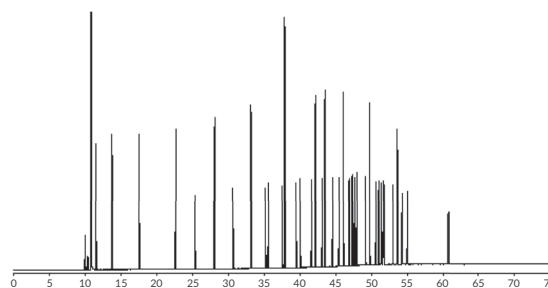
Fatty Ester Mixtures

Featured Product

Comprehensive Fatty Acid Methyl Ester Mixture 2

Item No. 29335

- Contains 37 saturated and mono- and polyunsaturated fatty acid methyl esters
- Intended for use as a reference standard in various sample types by GC-MS



Compare Fatty Ester Mixtures to Find the One Fit for Your Research

Learn more at www.caymanchem.com/FAME-mixtures

Fatty Amides

Unlabeled Standards

Item No.	Product Name
90350	Palmitoyl Ethanolamide
90050	Arachidonoyl Ethanolamide
90051	Arachidonoyl Glycine
90057	N-Arachidonoyl Dopamine
10005537	N-Arachidonoyl Taurine
10005455	N-Arachidonoyl-L-Serine
10964	Eicosapentaenoyl Ethanolamide

Popular matched pair items shown, over 165 fatty amides available online

Labeled Standards

Item No.	Product Name
10007824	Palmitoyl Ethanolamide-d ₄
10011178	Arachidonoyl Ethanolamide-d ₄
10007531	Arachidonoyl Glycine-d ₈
10007431	N-Arachidonoyl Dopamine-d ₈
26323	N-Arachidonoyl Taurine-d ₄
10007428	N-Arachidonoyl-L-Serine-d ₈
9001835	Eicosapentaenoyl Ethanolamide-d ₄

Fatty Acyl Carnitines

Unlabeled Standards

Item No.	Product Name
9001873	Propionyl-L-carnitine (chloride)
26555	Isovaleryl-L-carnitine (chloride)
26554	Hexanoyl-L-carnitine (chloride)
26558	Octanoyl-L-carnitine (chloride)
26550	Lauroyl-L-carnitine (chloride)
26556	Stearoyl-L-carnitine (chloride)
39014	Oleoyl-L-carnitine (chloride)

Popular matched pair items shown, over 55 fatty acyl carnitines available online

Labeled Standards

Item No.	Product Name
26579	Propionyl-L-carnitine-d ₃ (chloride)
26574	Isovaleryl-L-carnitine-d ₃ (chloride)
26570	Hexanoyl-L-carnitine-d ₃ (chloride)
26577	Octanoyl-L-carnitine-d ₃ (chloride)
26571	Lauroyl-L-carnitine-d ₃ (chloride)
26580	Stearoyl-L-carnitine-d ₃ (chloride)
26578	Oleoyl-L-carnitine-d ₃ (chloride)

Custom Synthesis & Mixtures

- Complex fatty acids
- Sphingolipids
- Stable isotope-labeled standards
- Glycerophospholipids
- Sterols
- Custom MaxSpec® mixtures

Learn more at www.caymanchem.com/chemicalsynthesis

Glycerolipids

Unlabeled Standards

Item No.	Product Name
23334	1,2,3-Tripalmitoyl Glycerol; TG(16:0/16:0/16:0)
16489	1,2,3-Triheptadecanoyl Glycerol; TG(17:0/17:0/17:0)
23335	1,2,3-Tristearoyl Glycerol; TG(18:0/18:0/18:0)
26871	1,2,3-Trioлеoyl Glycerol; TG(18:1/18:1/18:1)
10008650	1-Stearoyl-2-Arachidonoyl- <i>sn</i> -Glycerol; SAG
26827	1-Stearoyl- <i>rac</i> -glycerol; MG(18:0/0:0/0:0)
62160	2-Arachidonoyl Glycerol; 2-AG

Popular matched pair items shown, over 320 glycerolipids available online

Labeled Standards

Item No.	Product Name
28593	1,2,3-Tripalmitoyl-d ₃₁ Glycerol; TG(16:0/16:0/16:0)-d ₃₁
28918	1,2,3-Triheptadecanoyl Glycerol-d ₅ ; TG(17:0/17:0/17:0)-d ₅
28867	1,2,3-Tristearoyl Glycerol-d ₁₀₅ ; TG(18:0/18:0/18:0)-d ₁₀₅
41987	1,2,3-Trioлеoyl Glycerol-d ₅ ; TG(18:1/18:1/18:1)-d ₅
10009872	1-Stearoyl-2-Arachidonoyl-d ₈ - <i>sn</i> -Glycerol; SAG-d ₈
28919	1-Stearoyl- <i>rac</i> -glycerol-d ₄₀ ; MG(18:0-d ⁴⁰ /0:0/0:0)
362162	2-Arachidonoyl Glycerol-d ₅ ; 2-AG-d ₅

Glycerophospholipids

Unlabeled Standards

Item No.	Product Name
15097	1,2-Dimyristoyl- <i>sn</i> -glycero-3-PC; DMPC
10009473	1,2-Dipalmitoyl- <i>sn</i> -glycero-3-PC; DPPC
15092	1,2-Dipalmitoyl- <i>sn</i> -glycero-3-PE; DPPE
15086	1,2-Dipalmitoyl- <i>sn</i> -glycero-3-PG (sodium salt); DPPG
15100	1,2-Distearoyl- <i>sn</i> -glycero-3-PC; DSPC
25658	1-Palmitoyl-2-Arachidonoyl- <i>sn</i> -glycero-3-PC; PAPC
10010094	1-Palmitoyl-2-hydroxy- <i>sn</i> -glycero-3-PA; LPA 16:0
10172	1-Palmitoyl-2-hydroxy- <i>sn</i> -glycero-3-PC; LPC 16:0
10009864	1-Stearoyl-2-Arachidonoyl- <i>sn</i> -glycero-3-PC; PC(18:0/20:4)
25871	1-Stearoyl-2-Arachidonoyl- <i>sn</i> -glycero-3-PE; PE(18:0/20:4)
15107	1-Stearoyl-2-hydroxy- <i>sn</i> -glycero-3-PC; Lyso-PC
60900	PAF C-16
60906	Lyso-PAF C-16
60916	Lyso-PAF C-18

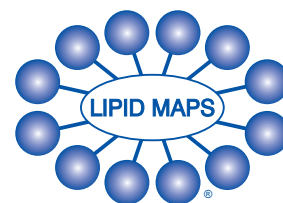
Popular matched pair items shown, over 280 glycerophospholipids available online

Labeled Standards

Item No.	Product Name
28753	1,2-Dimyristoyl-d ₅₄ - <i>sn</i> -glycero-3-PC; DMPC-d ₅₄
28154	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycero-3-PC; DPPC-d ₉
28155	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycero-3-PE; DPPE-d ₉
28953	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycero-3-PG; DPPG-d ₉
28752	1,2-Distearoyl-d ₇₀ - <i>sn</i> -glycero-3-PC; DSPC-d ₇₀
38384	1-Palmitoyl-d ₉ -2-Arachidonoyl- <i>sn</i> -glycero-3-PC; PAPC-d ₉
33479	1-Palmitoyl-d ₉ -2-hydroxy- <i>sn</i> -glycero-3-PA; LPA 16:0-d ₉
28153	1-Palmitoyl-d ₉ -2-hydroxy- <i>sn</i> -glycero-3-PC; LPC 16:0-d ₉
27928	1-Stearoyl-2-Arachidonoyl-d ₁₁ - <i>sn</i> -glycero-3-PC; PC(18:0/20:4-d ₁₁)
27929	1-Stearoyl-2-Arachidonoyl-d ₁₁ - <i>sn</i> -glycero-3-PE; PE(18:0/20:4-d ₁₁)
28230	1-Stearoyl-d ₃₅ -2-hydroxy- <i>sn</i> -glycero-3-PC; Lyso-PC-d ₃₅
360900	PAF C-16-d ₄
360906	Lyso-PAF C-16-d ₄
10010228	Lyso-PAF C-18 -d ₄

LIPID MAPS®

Cayman is a proud sponsor of the LIPID MAPS® consortium and supports its mission as a repository of data and protocols related to lipid research. Our catalog is increasingly aligned with the LIPID MAPS® classification and name recommendations, and we provide LIPID MAPS® with structures and information on new lipid standards available in our catalog.



Lipid Resource Center

Explore all products and resources for lipidomics and lipid analysis at www.caymanchem.com/lipids

Sphingolipids

Gangliosides

Unlabeled Standards

Item No.	Product Name
42979	C16 Ganglioside G _{M1} (d18:1/16:0) (ammonium salt)
42980	C18 Ganglioside G _{M1} (d18:1/18:0) (ammonium salt)
42981	C24:1 Ganglioside G _{M1} (d18:1/24:1) (ammonium salt)
15587	Ganglioside G _{M3} (bovine brain) (sodium salt)

Popular matched pair items shown, over 70 gangliosides available online

Glucosylceramides & Galactosylceramides

Unlabeled Standards

Item No.	Product Name
24352	C16 Lactosylceramide (d18:1/16:0)
40274	C16 α-Glucosylceramide (d18:1/16:0) (synthetic)
24863	C18 3'-sulfo Galactosylceramide (d18:1/18:0)
23210	C22 Glucosylceramide (d18:1/22:0)

Popular matched pair items shown, over 85 glucosylceramides and galactosylceramides available online

Globosides

Unlabeled Standards

Item No.	Product Name
24875	C16 Globotriaosylceramide (d18:1/16:0)
24877	C18 Globotriaosylceramide (d18:1/18:0)

Popular matched pair items shown, over 15 globosides available online

Ceramides, Sphingosines, & Sphingomyelins

Unlabeled Standards

Item No.	Product Name
62525	C6 Ceramide (d18:1/6:0)
10681	C16 Ceramide (d18:1/16:0)
19556	C18 Ceramide (d18:1/18:0)
62535	C24 Ceramide (d18:1/24:0)
62530	C24:1 Ceramide (d18:1/24:1(15Z))
10007907	Sphingosine (d18:1)
62570	Sphingosine-1-phosphate (d18:1)
10007946	C16 Sphingomyelin (d18:1/16:0)
20217	Phytosphingosine (yeast)

Popular matched pair items shown, over 225 ceramides, 180 sphingosines, and 45 sphingomyelins available online

Labeled Standards

Item No.	Product Name
31195	C16 Ganglioside G _{M1} -d ₉ (d18:1/16:0-d ₉) (ammonium salt)
24839	C18 Ganglioside G _{M1} -d ₃ (d18:1/18:0-d ₃) (ammonium salt)
42982	C24:1 Ganglioside G _{M1} -d ₁₈ (d18:1/24:1-d ₁₈) (ammonium salt)
39226	C18 Ganglioside G _{M3} -d ₃ (d18:1/18:0-d ₃) (ammonium salt, synthetic)

Labeled Standards

Item No.	Product Name
24625	C16 Lactosylceramide-d ₃ (d18:1/16:0-d ₃)
24621	C16 Glucosylceramide-d ₃ (d18:1/16:0-d ₃)
24624	C18 3'-sulfo Galactosylceramide-d ₃ (d18:1/18:0-d ₃)
9003462	C22 Glucosylceramide-d ₄ (d18:1/22:0-d ₄)

Labeled Standards

Item No.	Product Name
31199	C16 Globotriaosylceramide-d ₉ (d18:1/16:0-d ₉)
24626	C18 Globotriaosylceramide-d ₃ (d18:1/18:0-d ₃)

Labeled Standards

Item No.	Product Name
24826	C6 Ceramide-d ₇ (d18:1-d ₇ /6:0)
22787	C16 Ceramide-d ₇ (d18:1-d ₇ /16:0)
22788	C18 Ceramide-d ₇ (d18:1-d ₇ /18:0)
22789	C24 Ceramide-d ₇ (d18:1-d ₇ /24:0)
22790	C24:1 Ceramide-d ₇ (d18:1-d ₇ /24:1(15Z))
22786	Sphingosine-d ₇ (d18:1)
10008121	Sphingosine-1-phosphate-d ₇ (d18:1)
27551	C16 Sphingomyelin-d ₉ (d18:1/16:0-d ₉)
9004341	Phytosphingosine-d ₇

View Tips on Sphingolipid Solubilization, Isolation, Analysis, and Storage

Learn more from our lipid experts at [z](#)

Sterol Lipids

Cholesterol Esters

Unlabeled Standards

Item No.	Product Name
22595	Cholesteryl Arachidonate
22597	Cholesteryl Linoleate
26473	Cholesteryl Palmitate

Popular matched pair items shown, over 45 cholesteryl esters available online

Oxysterols

Unlabeled Standards

Item No.	Product Name
19518	4 β -hydroxy Cholesterol
25538	5 α ,6 β -Dihydroxycholestanol
20098	7 α -hydroxy Cholesterol
25442	7 α -hydroxy-4-Cholesten-3-one
20099	7 β -hydroxy Cholesterol
11097	25-hydroxy Cholesterol

Popular matched pair items shown, over 15 oxysterols available online

Labeled Standards

Item No.	Product Name
39085	Cholesteryl Arachidonate-d ₈
37920	Cholesteryl Linoleate-d ₁₁
28123	Cholesteryl Palmitate-d ₉

Labeled Standards

Item No.	Product Name
25971	4 β -hydroxy Cholesterol-d ₇
25968	5 α ,6 β -Dihydroxycholestanol-d ₇
25547	7 α -hydroxy Cholesterol-d ₇
26100	7 α -hydroxy-4-Cholesten-3-one-d ₇
25548	7 β -hydroxy Cholesterol-d ₇
11099	25-hydroxy Cholesterol-d ₆

Oxysterol Derivatization MaxSpec® Kit - Item No. 601540

An oxidation and derivatization kit for mass spec analysis of oxysterols

Bile Acids

Unlabeled Standards

Item No.	Product Name
20250	Cholic Acid; CA
16215	Taurocholic Acid (sodium salt); TCA
20276	Glycocholic Acid; GCA
10011286	Chenodeoxycholic Acid; CDCA
20275	Taurochenodeoxycholic Acid (sodium salt); TCDCA
16942	Glycochenodeoxycholic Acid (sodium salt hydrate); GCDCA
20253	Lithocholic Acid; LCA
20756	Deoxycholic Acid; DCA
15935	Taurodeoxycholic Acid (sodium salt hydrate); TDCA

Popular matched pair items shown, over 155 bile acids available online

Labeled Standards

Item No.	Product Name
20849	Cholic Acid-d ₄ ; CA-d ₄
21891	Taurocholic Acid-d ₄ (sodium salt); TCA-d ₄
21889	Glycocholic Acid-d ₄ ; GCA-d ₄
20848	Chenodeoxycholic Acid-d ₄ ; CDCA-d ₄
21905	Taurochenodeoxycholic Acid-d ₄ (sodium salt); TCDCA-d ₄
21890	Glycochenodeoxycholic Acid-d ₄ ; GCDCA-d ₄
20831	Lithocholic Acid-d ₄ ; LCA-d ₄
20851	Deoxycholic Acid-d ₄ ; DCA-d ₄
32983	Taurodeoxycholic Acid-d ₄ (sodium salt); TDCA-d ₄

MaxSpec® Mixtures for Bile Acids

Contains primary and secondary bile acids, as well as glycine- and taurine-conjugated bile acids

Item No.	Product Name
33505	Bile Acids MaxSpec® Discovery Mixture

Item No.	Product Name
33506	Deuterated Bile Acids MaxSpec® Discovery Mixture