

MaxSpec® Mixtures for LC-MS

Cayman's MaxSpec® mixtures streamline workflows by reducing the effort and variability of mixing individual internal and calibration standards. Each mixture undergoes batch-specific LC-MS content analysis and ongoing stability testing to ensure quality.



MaxSpec® Quality Guarantee

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



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

Bile Acids

Item No.	Product Name	Contains (1,200-4,300 nM of each compound)
33505	Bile Acids MaxSpec® Discovery Mixture*	Tauroursodeoxycholic Acid, Taurocholic Acid, Taurochenodeoxycholic Acid, Taurodeoxycholic Acid, Taurothiocholic Acid, Chenodeoxycholic Acid, Hyodeoxycholic Acid, Lithocholic Acid, Deoxycholic Acid, Glycocholic Acid, Glycoursodeoxycholic Acid, Ursodeoxycholic Acid, Glycochenodeoxycholic Acid, Glycolithocholic Acid, Glycodeoxycholic Acid, Glycohyodeoxycholic Acid, Taurohyodeoxycholic Acid, Cholic Acid

*Also available: Deuterated Bile Acids MaxSpec® Discovery Mixture (Item No. 33506)

Bioactive Eicosanoids, Precursors, & Metabolites

Item No.	Product Name	Contains (1 µg/ml of each compound)
19101	Primary COX and LOX MaxSpec® LC-MS Mixture*	6-keto Prostaglandin F _{1α} , Thromboxane B ₂ , Prostaglandin F _{2α} , Prostaglandin E ₂ , Prostaglandin D ₂ , 12(S)-HHTrE, 15(S)-HETE, 12(S)-HETE, 5(S)-HETE
19300	Primary Prostaglandin Metabolite MaxSpec® LC-MS Mixture*	11β-Prostaglandin F _{2α} , 15-keto Prostaglandin F _{2α} , 15-keto Prostaglandin E ₂ , 13,14-dihydro-15-keto Prostaglandin F _{2α} , 13,14-dihydro-15-keto Prostaglandin E ₂ , 13,14-dihydro-15-keto Prostaglandin D ₂ , Prostaglandin J ₂ , 15-deoxy-Δ ^{12,14} -Prostaglandin J ₂
19422	Secondary Prostaglandin Metabolite MaxSpec® LC-MS Mixture	tetranor-PGFM, tetranor-PGEM, tetranor-PGDM, 13,14-dihydro-15-keto Prostaglandin F _{2α} , 13,14-dihydro-15-keto Prostaglandin E ₂ , 13,14-dihydro-15-keto Prostaglandin D ₂

*Also available: Deuterated Primary COX and LOX MaxSpec® LC-MS Mixture (Item No. 19228) and Deuterated Primary Prostaglandin Metabolite MaxSpec® LC-MS Mixture (Item No. 19304)

Custom Formulations and Mixtures Available
Email sales@caymanchem.com to discuss your specific project needs

Dietary & Microbiome-Related Fatty Acids

Item No.	Product Name	Contains (10 µg/ml of each compound)
18824	Short-chain Fatty Acid MaxSpec® LC-MS Mixture	Propionic Acid (C3:0), Butyric Acid (C4:0), Hexanoic Acid (C6:0), Heptanoic Acid (C7:0), Octanoic Acid (C8:0), Nonanoic Acid (C9:0), Decanoic Acid (C10:0), Dodecanoic Acid (C12:0)
19227	Odd-chain Fatty Acid MaxSpec® LC-MS Mixture	Undecanoic Acid (C11:0), Tridecanoic Acid (C13:0), 14-Pentadecenoic Acid (C15:1), Pentadecanoic Acid (C15:0), <i>cis</i> -10-Heptadecenoic Acid (C17:1), Heptadecanoic Acid (C17:0), <i>cis</i> -10-Nonadecenoic Acid (C19:1), Nonadecanoic Acid (C19:0)
17941	Polyunsaturated Fatty Acid MaxSpec® LC-MS Mixture	Adrenic Acid (C22:4), Arachidonic Acid (C20:4), Dihomo-γ-Linolenic Acid (C20:3), Docosahexaenoic Acid (C22:6), Docosapentaenoic Acid (C22:5), Eicosapentaenoic Acid (C20:5), Linoleic Acid (C18:2), α-Linolenic Acid (C18:3 n-3), γ-Linolenic Acid (C18:3 n-6), Stearidonic Acid (C18:4)
17942	Saturated/Monounsaturated Fatty Acid MaxSpec® LC-MS Mixture	Arachidic Acid (C20:0), Lauric Acid (C12:0), Lignoceric Acid (C24:0), Myristic Acid (C14:0), Nervonic Acid (C24:1), Oleic Acid (C18:1), Palmitic Acid (C16:0), Palmitoleic Acid (C16:1), Stearic Acid (C18:0)

Oxylipins

Item No.	Product Name	Contains (1-10 µg/ml of each compound)
22638	ALA and GLA Oxylipin MaxSpec® LC-MS Mixture	9(S)-HOTrE, 13(S)-HOTrE, 13(S)-HOTrE(γ)
20666	Arachidonic Acid Oxylipin MaxSpec® LC-MS Mixture*	15(S)-HETE, 11(S)-HETE, 15-OxoETE, 8(S)-HETE, 12(S)-HETE, 9(R)-HETE, 5(S)-HETE, 12-OxoETE, 5-OxoETE
22280	DHA Oxylipin MaxSpec® LC-MS Mixture	(±)20-HDHA, (±)16-HDHA, 17(S)-HDHA, (±)13-HDHA, (±)10-HDHA, 14(S)-HDHA, (±)11-HDHA, (±)7-HDHA, (±)8-HDHA, (±)4-HDHA
22281	Dihomo-γ-Linolenic Acid Metabolite MaxSpec® LC-MS Mixture	Prostaglandin F _{1α} , Prostaglandin E ₁ , 15-keto Prostaglandin E ₁ , 13,14-dihydro-15-keto Prostaglandin F _{1α} , 13,14-dihydro-15-keto Prostaglandin E ₁ , 8(S)-HETrE, 15(S)-HETrE, 12(S)-HETrE
21393	EPA Oxylipin MaxSpec® LC-MS Mixture	Prostaglandin E ₃ , Prostaglandin F _{3α} , (±)18-HEPE, 15(S)-HEPE, (±)8-HEPE, (±)11-HEPE, (±)9-HEPE, 12(S)-HEPE, 5(S)-HEPE
22640	Leukotriene B ₄ Pathway MaxSpec® LC-MS Mixture	18-carboxy dinor Leukotriene B ₄ , 20-carboxy Leukotriene B ₄ , 20-hydroxy Leukotriene B ₄ , Leukotriene B ₄ , 5(S)-HETE
20794	Linoleic Acid Oxylipins MaxSpec® LC-MS Mixture*	(±)12(13)-DiHOME, (±)9(10)DiHOME, 9(S)-HODE, 13(S)-HODE, 13-OxoODE, 9-OxoODE, (±)12(13)-EpOME, (±)9(10)-EpOME

*Also available: Deuterated Arachidonic Acid Oxylipin MaxSpec® LC-MS Mixture (Item No. 20668) and Deuterated Linoleic Acid Oxylipins MaxSpec® LC-MS Mixture (Item No. 20795)

CYP450 Metabolites

Item No.	Product Name	Contains (1-10 µg/ml of each compound)
20665	Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture*	(±)8(9)-EET, (±)11(12)-EET, (±)14(15)-EET, (±)5(6)-DiHET, (±)8(9)-DiHET, (±)11(12)-DiHET, (±)14(15)-DiHET, 19(R)-HETE, 20-HETE
22639	Docosahexaenoic Acid CYP450 Oxylipins MaxSpec® LC-MS Mixture	(±)19(20)-DiHDPA, (±)16(17)-DiHDPA, (±)13(14)-DiHDPA, (±)10(11)-DiHDPA, (±)7(8)-DiHDPA, (±)19(20)-EpDPA, (±)16(17)-EpDPA, (±)13(14)-EpDPA, (±)10(11)-EpDPA, (±)7(8)-EpDPA
21394	EPA CYP450 Oxylipin MaxSpec® LC-MS Mixture	(±)17(18)-DiHETE, (±)11(12)-DiHETE, (±)14(15)-DiHETE, (±)8(9)-DiHETE, (±)5(6)-DiHETE, (±)17(18)-EpETE, (±)14(15)-EpETE, (±)8(9)-EpETE, (±)11(12)-EpETE

*Also available: Deuterated Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture (Item No. 20667)

Specialized Pro-Resolving Mediators

Item No.	Product Name	Contains (100 ng/ml of each compound)
19412	Lipoxin MaxSpec® LC-MS Mixture	Lipoxin B ₄ , Lipoxin A ₄ , 15(R)-Lipoxin A ₄ , Arachidonic Acid
18702	SPM D-series MaxSpec® LC-MS Mixture	Resolvin D3, Resolvin D2, 17(R)-Resolvin D1, Resolvin D1, Resolvin D5, Docosahexaenoic Acid
19417	SPM E-series MaxSpec® LC-MS Mixture	Resolvin E1, (±)18-HEPE, Eicosapentaenoic Acid

Vasoactive Eicosanoids

Item No.	Product Name	Contains (1 µg/ml of each compound)
19667	Primary Vascular Eicosanoid MaxSpec® LC-MS Mixture	6-keto Prostaglandin F _{1α} , Thromboxane B ₂ , 11-dehydro Thromboxane B ₂ , 12(S)-HHTrE, 20-HETE, 12(S)-HETE, 12(S)-HETrE, (±)14(15)-EET, (±)11(12)-EET, (±)8(9)-EET

Also available: Vascular Eicosanoid Urinary Metabolite MaxSpec® LC-MS Mixture (Item No. 19668)

Oxidative Stress

Item No.	Product Name	Contains (1 µg/ml of each compound)
18701	Oxidative Stress MaxSpec® LC-MS Mixture	8-Hydroxyguanosine, 8-Hydroxy-2'-deoxyguanosine, 2,3-dinor-8-iso Prostaglandin F _{2α} , 8-iso Prostaglandin F _{1α} , 8-iso Prostaglandin F _{2α} , (±)5-iPF _{2α} -VI, 8-iso-15-keto Prostaglandin F _{2α} , 8-iso Prostaglandin E ₂ , 8,12-iso-iPF _{2α} -VI