

PRODUCT INFORMATION



EPA Oxylin MaxSpec® LC-MS Mixture Item No. 21393

Supplied as: A solution in ethanol (1 µg/ml each compound)
Fill volume: >1 ml
Storage: -20°C
Stability: ≥5 years

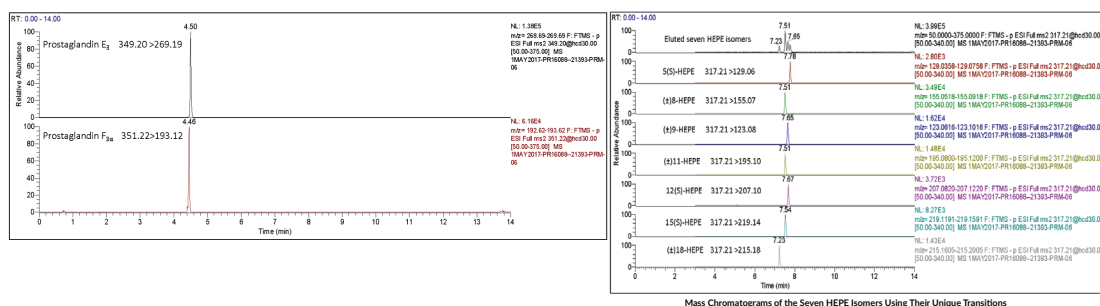
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

EPA oxylin MaxSpec® LC-MS mixture contains oxylin metabolites derived from eicosapentaenoic acid (EPA; Item Nos. 90110 | 21908). The mixture is supplied in an amber ampule in which the headspace has been purged with argon to prevent lipid oxidation. This product has been designed for direct use in LC-MS applications. The solution may be used as a system suitability standard or tuning standard. After opening, we recommend that the mixture be transferred immediately to a 1 ml glass screw cap vial, to prevent solvent evaporation, and stored at -20°C. The mixture should be discarded after multiple freeze/thaw cycles.

This mixture contains Prostaglandin E₃ (Item No. 14990), Prostaglandin F_{3α} (Item No. 16990), 5(S)-HEPE (Item No. 32210), (±)8-HEPE (Item No. 32340), (±)9-HEPE (Item No. 32400), (±)11-HEPE (Item No. 32500), 12(S)-HEPE (Item No. 32550), 15(S)-HEPE (Item No. 32710), and (±)18-HEPE (Item No. 32840).

Contents



Item Number: 21393		EPA Oxylin MaxSpec® LC-MS Mixture			
Item Number	Item Name	Formula	Mass (Da)	Transitions (m/z):	RT (min)
14990	Prostaglandin E ₃	C ₂₀ H ₃₀ O ₅	350.5	349.20>269.19	4.50
16990	Prostaglandin F _{3α}	C ₂₀ H ₃₂ O ₅	352.5	351.22>193.12	4.46
32210	5(S)-HEPE	C ₂₀ H ₃₀ O ₃	318.5	317.21>129.06	7.78
32340	(±)8-HEPE	C ₂₀ H ₃₀ O ₃	318.5	317.21>155.07	7.51
32400	(±)9-HEPE	C ₂₀ H ₃₀ O ₃	318.5	317.21>123.08	7.65
32500	(±)11-HEPE	C ₂₀ H ₃₀ O ₃	318.5	317.21>195.10	7.51
32550	12(S)-HEPE	C ₂₀ H ₃₀ O ₃	318.5	317.21>207.10	7.67
32710	15(S)-HEPE	C ₂₀ H ₃₀ O ₃	318.5	317.21>219.14	7.54
32840	(±)18-HEPE	C ₂₀ H ₃₀ O ₃	318.5	317.21>215.18	7.23
LC-MS/MS: Ultimate 3000 UHPLC&Q-Exactive Quadrupole-Orbitrap (Thermo Scientific)					
Mobile Phase A: Water + 0.1% Formic Acid					
Mobile Phase B: Acetonitrile + 0.1% Formic Acid					
Column: Waters BEH C8, 2.1 x 100 mm, 1.7 µm			Flow Rate: 400 µl/min		
Negative Electrospray Ionization			Parallel Reaction Monitoring (PRM)		

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/14/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM