

PRODUCT INFORMATION



EPA CYP450 Oxylipin MaxSpec® LC-MS Mixture Item No. 21394

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

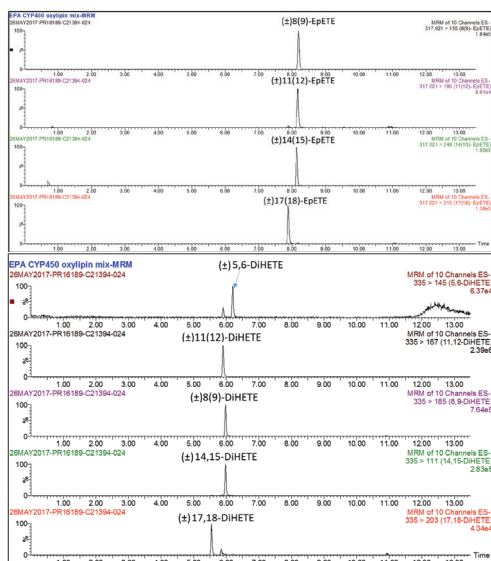
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

The EPA CYP450 oxylipin MaxSpec® LC-MS mixture contains various regioisomers of epoxides and dihydroxy fatty acids derived from the sequential metabolism of eicosapentaenoic acid (EPA; Item Nos. 90110 | 21908) via CYP450 and soluble epoxide hydrolase. 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 (±)8(9)-EpETE (Item No. 10470), (±)11(12)-EpETE (Item No. 10462), (±)14(15)-EpETE (Item No. 10173), (±)17(18)-EpETE (Item No. 50861), (±)5(6)-DiHETE (Item No. 10467), (±)11(12)-DiHETE (Item No. 10466), (±)8(9)-DiHETE (Item No. 10473), (±)14(15)-DiHETE (Item No. 10006998), and (±)17(18)-DiHETE (Item No. 10006999).

Contents



Item Number: 21394		EPA CYP450 Oxylipin MaxSpec® LC-MS Mixture			
Item Number	Item Name	Formula	Mass (Da)	Transition (m/z)	RT (min)
10470	(±)8(9)-EpETE	C ₂₀ H ₃₀ O ₃	318.5	317>155	8.2
10462	(±)11(12)-EpETE	C ₂₀ H ₃₀ O ₃	318.5	317>195	8.2
10173	(±)14(15)-EpETE	C ₂₀ H ₃₀ O ₃	318.5	317>248	8.1
50861	(±)17(18)-EpETE	C ₂₀ H ₃₀ O ₃	318.5	317>215	7.9
10467	(±)5(6)-DiHETE	C ₂₀ H ₃₂ O ₄	336.5	335>145	6.2
10466	(±)11(12)-DiHETE	C ₂₀ H ₃₂ O ₄	336.5	335>167	5.9
10473	(±)8(9)-DiHETE	C ₂₀ H ₃₂ O ₄	336.5	335>185	6.0
10006998	(±)14(15)-DiHETE	C ₂₀ H ₃₂ O ₄	336.5	335>111	5.9
10006999	(±)17(18)-DiHETE	C ₂₀ H ₃₂ O ₄	336.5	335>203	5.5
LC-MS: Waters Acquity UPLC-Xevo TQ-S Micro					
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: 300 µl/min		
Negative Electrospray Ionization			MRM Scan		

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