

PRODUCT INFORMATION



Arachidonic Acid Oxylin LC-MS Mixture

Item No. 20666

Supplied as: A solution in ethanol (10 µg/ml each compound)

Storage: -80°C

Stability: ≥3 years

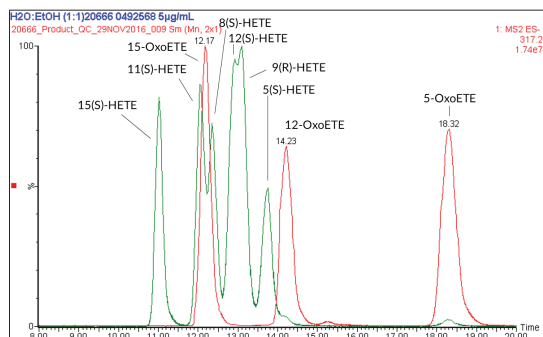
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

The arachidonic acid oxylin LC-MS mixture contains oxylin metabolites derived from arachidonic acid (Item No 90010). 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 serially diluted for preparation of calibrators and QC standards and/or used directly 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 5(S)-HETE (Item No. 34230), 5-OxoETE (Item No. 34250), 8(S)-HETE (Item No. 34360), 9(R)-HETE (Item No. 34405), 11(S)-HETE (Item No. 34510), 12(S)-HETE (Item No. 34570), 12-OxoETE (Item No. 34580), 15(S)-HETE (Item No. 34720), and 15-OxoETE (Item No. 34730).

Contents



Item Number: 20666		Arachidonic Acid Oxylin LC-MS Mixture		
Item Number	Item Name	Formula Weight	MS/MS Transition	RT (min)
34720	15(S)-HETE	320.5	319>219	11.08
34510	11(S)-HETE	320.5	319>195	12.03
34730	15-OxoETE	318.5	317>113	12.17
34360	8(S)-HETE	320.5	319>155	12.42
34570	12(S)-HETE	320.5	319>179	12.87
34405	9(R)-HETE	320.5	319>139	13.08
34230	5(S)-HETE	320.5	319>115	13.77
34580	12-OxoETE	318.5	317>153	14.23
34250	5-OxoETE	318.5	317>203	18.32

LC-MS Conditions:

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
LC Gradient: Isocratic 55%B over 20 minutes	
Negative Electrospray Ionization	

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, 12/18/2018

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