

PRODUCT INFORMATION



Deuterated Arachidonic Acid Oxylipin MaxSpec® LC-MS Mixture Item No. 20668

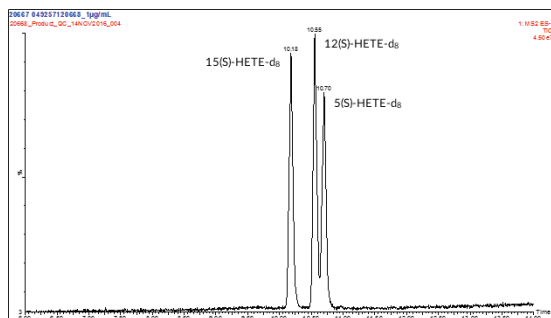
Supplied as: A solution in acetonitrile (1 µg/ml each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥5 years

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

Description and Contents

The deuterated arachidonic acid oxylipin MaxSpec® LC-MS Mixture contains deuterated forms of oxylipin 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. This mixture is ideally suited for use as an internal standard in mass spectrometry applications. 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-d₈ (Item No. 334230), 12(S)-HETE-d₈ (Item No. 334570), and 15(S)-HETE-d₈ (Item No. 334720). The unmodified forms are involved in the immune response.



Item Number: 20668		Deuterated Arachidonic Acid Oxylipin MaxSpec® LC-MS Mixture	
Item Number	Item Name	Formula Weight:	MS/MS Transition:
334720	15(S)-HETE-d ₈	328.5	327>226
334570	12(S)-HETE-d ₈	328.5	327>184
334230	5(S)-HETE-d ₈	328.5	327>116
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: 25% B to 50% B over 7.5 minutes, 50% to 60% 12.5 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, 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