

PRODUCT INFORMATION



Deuterated Linoleic Acid Oxylipins MaxSpec[®] LC-MS Mixture Item No. 20795

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

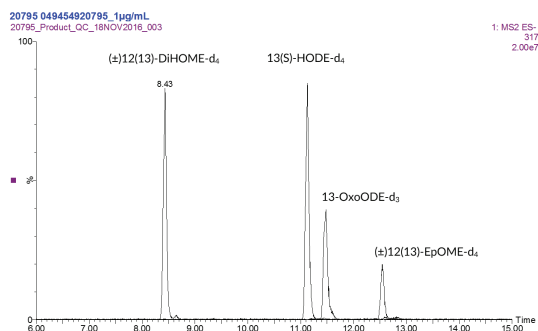
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

This mixture contains deuterated versions of oxylipin metabolites derived from linoleic acid (Item No. 90150). 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 deuterated forms of (±)12(13)-DiHOME (Item No. 10009832), 13(S)-HODE (Item No. 38610), 13-OxoODE (Item No. 38620), and (±)12(13)-EpOME (Item No. 52450), each of which have been shown to be involved in the physiological or pathological responses related to cell proliferation or oxidative burst and are potential markers for (or potential contributors to) various human diseases.

Contents



Item Number: 20795		Deuterated Linoleic Acid Oxylipins MaxSpec [®] LC-MS Mixture		
Item Number	Item Name	Formula Weight:	MS/MS Transition:	RT (min)
10009994	(±)12(13)-DiHOME-d ₄	318.5	317>185	8.43
338610	13(S)-HODE-d ₄	300.5	299>198	11.13
338620	13-OxoODE-d ₃	297.5	296>198	11.43
10009996	(±)12(13)-EpOME-d ₄	300.5	299>198	12.55

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, 25°C	Flow Rate: 400 µl/min
LC Gradient: Isocratic 25%B to 95%B over 24 minutes	
Negative Electrospray Ionization	Full MS/Survey 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, 02/08/2024

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