

PRODUCT INFORMATION



DHA Oxylin MaxSpec® LC-MS Mixture

Item No. 22280

Supplied as: A solution in ethanol (1 µg/ml each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥7 years

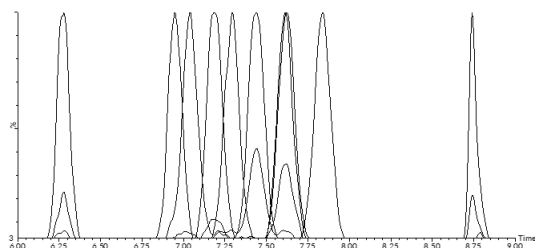
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

The DHA Oxylin MaxSpec® LC-MS Mixture contains oxylin metabolites derived from docosahexaenoic acid (DHA; Item No. 90310). 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 (±)4-HDHA (Item No. 33200), (±)7-HDHA (Item No. 33300), (±)8-HDHA (Item No. 33350), (±)10-HDHA (Item No. 33400), (±)11-HDHA (Item No. 33450), (±)13-HDHA (Item No. 33500), 14(S)-HDHA (Item No. 15253), (±)16-HDHA (Item No. 33600), 17(S)-HDHA (Item No. 10009799), and (±)20-HDHA (Item No. 33750).

Contents



Item Number: 22280		DHA Oxylin MaxSpec® LC-MS Mixture		
Item Number	Item Name	Formula Weight	MS/MS Transition	RT (min)
33200	(±)4-HDHA	344.5	343>101	8.74
33300	(±)7-HDHA	344.5	343>141	7.62
33350	(±)8-HDHA	344.5	343>125	7.84
33400	(±)10-HDHA	344.5	343>181	7.29
33450	(±)11-HDHA	344.5	343>165	7.61
33500	(±)13-HDHA	344.5	343>221	7.19
15253	14(S)-HDHA	344.5	343>205	7.44
33600	(±)16-HDHA	344.5	343>261	6.94
10009799	17(S)-HDHA	344.5	343>245	7.04
33750	(±)20-HDHA	344.5	343>285	6.28

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 10%B to 50%B over 1 minute, 7 minute hold, 95% B wash	
Negative Electrospray Ionization	MRM

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, 07/17/2023

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