

# PRODUCT INFORMATION



## Docosahexaenoic Acid CYP450 Oxylinps LC-MS Mixture

Item No. 22639

**Supplied as:** A solution in ethanol (1 µg/ml of each compound)

**Storage:** -20°C

**Stability:** ≥2 years

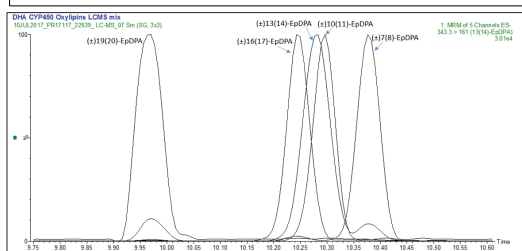
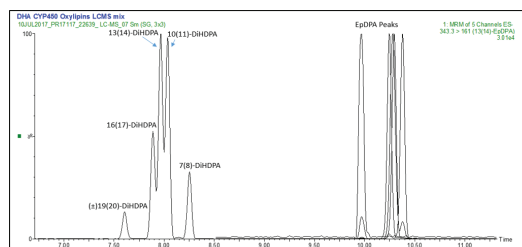
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

The docosahexaenoic acid (DHA) CYP450 oxylinps LC-MS mixture contains oxylinp metabolites derived from 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 (±)7(8)-EpDPA (Item No. 10465), (±)10(11)-EpDPA (Item No. 10471), (±)13(14)-EpDPA (Item No. 10464), (±)16(17)-EpDPA (Item No. 10174), (±)19(20)-EpDPA (Item No. 10175), 7(8)-DiHDPA (Item No. 10469), 10(11)-DiHDPA (Item No. 18174), 13(14)-DiHDPA (Item No. 18175), 16(17)-DiHDPA (Item No. 18176), and (±)19(20)-DiHDPA (Item No. 10007001).

### Contents



Item Number: 22639		DHA CYP450 Oxylinps LC-MS Mixture		
Item Number	Item Name	Mass (Da)	M/Z Transition	RT (min)
10007001	(±)19(20)-DiHDPA	362.5	361.5>229.4	7.61
18176	16(17)-DiHDPA	362.5	361.5>233.1	7.89
18175	13(14)-DiHDPA	362.5	361.3>193.1	7.97
18174	10(11)-DiHDPA	362.5	361.3>153.2	8.01
10469	7(8)-DiHDPA	362.5	361.3>127.1	8.25
10175	(±)19(20)-EpDPA	344.5	343.5>241.6	9.98
10174	(±)16(17)-EpDPA	344.5	343.3>274.4	10.23
10464	(±)13(14)-EpDPA	344.5	343.3>161.3	10.27
10471	(±)10(11)-EpDPA	344.5	343.4>153.2	10.30
10465	(±)7(8)-EpDPA	344.5	343.5>112.8	10.37

LC-MS/MS: Waters Acquity UPLC-XEVO TQD  
 Mobile Phase: A: Water + 0.1% Formic Acid; B: Acetonitrile + 0.1% Formic Acid

Column: Waters BEH C8, 2.1 x 100 mm, 1.7 µm. Negative Electrospray Ionization	Flow Rate: 400 µl/min 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, 03/12/2019

#### 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