

PRODUCT INFORMATION



Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture Item No. 20665

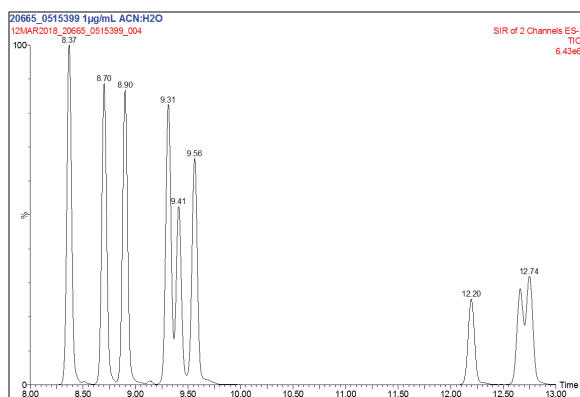
Supplied as: A solution in ethanol (10 µg/ml each compound)
Fill Volume: >1 ml
Storage: -20°C
Stability: ≥5 years

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

Description

The Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture contains CYP450 metabolites derived from arachidonic acid. 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.

Contents



Item Number: 20665			Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture	
Item Number	Item Name	Retention Time (min)	Formula Weight	MS/MS Transition
51651	(±)14(15)-DIHET	8.37	338.5	337>207
51511	(±)11(12)-DIHET	8.70	338.5	337>167
51351	(±)8(9)-DIHET	8.90	338.5	337>185
51211	(±)5(6)-DIHET	9.31	338.5	337>145
10007767	19(R)-HETE	9.41	320.5	319>231
90030	20-HETE	9.56	320.5	319>245
50651	(±)14(15)-EET	12.20	320.5	319>219
50511	(±)11(12)-EET	12.66	320.5	319>167
50351	(±)8(9)-EET	12.74	320.5	319>127
LC-MS Conditions: Waters Acquity UPLC, Xevo TQD				
Mobile Phase A: Water + 0.1% Formic acid		Time (min)	%A	%B
Mobile Phase B: Acetonitrile + 0.1% Formic acid		0.00	75	25
Column: Waters BEH C8, 2.1 x 100 mm, 1.7 µm		7.50	50	50
Flow Rate: 400 µl/min		12.50	40	60
Negative Electrospray Ionization		15.00	40	60
SIR Scan		15.01	5	95
		16.00	5	95
		16.01	75	25

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