

# PRODUCT INFORMATION

## Deuterated Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture

Item No. 20667

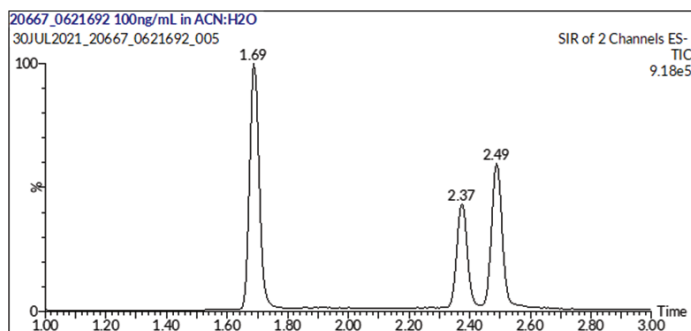
**Supplied as:** A solution in ethanol (1 µg/ml each compound)  
**Fill Volume:** >1 ml  
**Storage:** -20°C  
**Stability:** ≥5 years

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

### Description and Contents

The deuterated arachidonic acid CYP450 metabolite LC-MS mixture contains deuterated versions of HETE and EET metabolites produced by the cytochrome P450 pathway of arachidonic acid metabolism. 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 20-HETE (Item No. 90030), (±)11(12)-EET (Item No. 50511), and (±)14(15)-EET (Item No. 50651), each of which have been shown to be involved in either vasoconstriction or hypertension.



Item Number: 20667 Deuterated Arachidonic Acid CYP450 Metabolite MaxSpec® LC-MS Mixture			
Item No.	Item Name	Formula Weight	Retention Time (min)
390030	20-HETE-d <sub>8</sub>	326.5	1.69
10006413	(±)14(15)-EET-d <sub>11</sub>	331.5	2.37
10006410	(±)11(12)-EET-d <sub>11</sub>	331.5	2.49
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: 15%B to 50%B over 12 min, 50%B to 95%B over 1 min			
Negative Electrospray Ionization			SIR 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, 04/04/2024

#### 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