

# PRODUCT INFORMATION

## Oxylipins MaxSpec® Discovery Mixture 2 at 5 µg/ml each

Item No. 33597

**Supplied as:** A mixture in 1 ml of ethanol:methyl acetate (75:25); in a deactivated glass ampule  
**Concentration:** See certificate of analysis for verified concentrations  
**Storage:** -20°C  
**Stability:** ≥3 years; *Stability testing is ongoing to ensure concentration accuracy. The certificate of analysis and product expiry date will be updated upon completion of testing.*  
**Special Conditions:** Store upright and unopened at -20°C. Warm to room temperature prior to opening. Light sensitive.

### Description

Oxylipins MaxSpec® Discovery Mixture 2 contains 10 oxylipins derived from cytochrome P450-mediated metabolism of either arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607) or linoleic acid (Item Nos. 90150 | 90150.1 | 21909). It also contains a dehydrogenase metabolite of 5(S)-HETE (Item No. 34230) and two metabolites produced by the non-enzymatic hydrolysis of leukotriene A<sub>4</sub>.

Oxylipins MaxSpec® Discovery Mixture 2 has been prepared specifically for mass spectrometry or any application where quantitative reproducibility is required. The solution has been prepared gravimetrically and is supplied in a deactivated glass ampule sealed under argon. The concentration was verified by comparison to an independently prepared calibration standard. The verified concentration is provided on the certificate of analysis. This Oxylipins MaxSpec® Discovery Mixture 2 is guaranteed to meet identity, purity, stability, and concentration specifications and is provided with a batch-specific certificate of analysis. Ongoing stability testing is performed to ensure the concentration remains accurate throughout the shelf life of the product.

**Note:** The amount of solution added to the vial is in excess of the listed amount. Therefore, it is necessary to accurately measure volumes for preparation of calibration standards. Follow recommended storage and handling conditions to maintain product quality.

### Contents

The Oxylipins MaxSpec® Discovery Mixture 2 contains the following:

Item No.	Item Name (Concentration: 5 µg/ml each)
31028	(±)12(13)-EpOME MaxSpec® Standard
31027	(±)12(13)-DiHOME MaxSpec® Standard
31026	(±)9(10)-EpOME MaxSpec® Standard
31025	(±)9(10)-DiHOME MaxSpec® Standard
10007253	5(S),6(S)-DiHETE MaxSpec® Standard
10007252	5(S),6(R)-DiHETE MaxSpec® Standard
10007244	5-OxoETE MaxSpec® Standard
10007267	(±)14(15)-DiHET MaxSpec® Standard
10007265	(±)8(9)-DiHET MaxSpec® Standard
10007264	(±)5(6)-DiHET MaxSpec® Standard
10007263	(±)14(15)-EET MaxSpec® Standard
10007262	(±)11(12)-EET MaxSpec® Standard
10007261	(±)8(9)-EET MaxSpec® Standard

**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, 03/21/2023

**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