

PRODUCT INFORMATION

Deuterated Lipidomics MaxSpec® Mixture

Item No. 40974

Supplied as: A mixture in chloroform:methanol:ethanol (6:4:1); in a deactivated glass ampule
Fill Volume: >1 ml
Concentration: See certificate of analysis for verified concentrations
Storage: -80°C
Stability: ≥1 year; *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 -80°C. Warm to room temperature prior to opening. Light sensitive.

Description

Deuterated Lipidomics MaxSpec® Mixture (Item No. 40974) is a MaxSpec® mixture containing deuterated versions of various lipid species. The mixture contains deuterated glycerolipid, sphingolipid, sterol lipid, fatty acid, fatty ester, and glycerophospholipid species. The mixture is supplied in an amber ampule in which the headspace has been purged with argon. It is designed for use in method development and for direct analysis of these lipid species using LC-MS and LC-MS/MS instrumentation. It is ideally suited as a multiplexed internal standard for lipidomics analysis.

Deuterated Lipidomics MaxSpec® Mixture has been prepared specifically for mass spectrometry and related applications 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 Deuterated Lipidomics MaxSpec® Mixture 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.*

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, 05/20/2025

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

PRODUCT INFORMATION



Contents

The Deuterated Lipidomics MaxSpec® Mixture contains the following:

Item No.	Item Name	Concentration (µg/ml)
27588	1-Palmitoyl-d ₉ -2-hydroxy- <i>sn</i> -glycero-3-PE	10
28153	1-Palmitoyl-d ₉ -2-hydroxy- <i>sn</i> -glycero-3-PC	10
9003263	Heptadecanoyl-L-carnitine-d ₃ (chloride)	5
30557	Palmitic Acid-d ₅	40
27551	C16 Sphingomyelin-d ₉ (d18:1/16:0-d ₉)	10
25265	Cholesterol-d ₇	100
28156	PtdIns-(1-palmitoyl-d ₉ , 2-palmitoyl) (sodium salt)	10
28152	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycero-3-PS	10
28953	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycero-3-PG	10
22787	C16 Ceramide-d ₇ (d18:1-d ₇ /16:0)	5
28155	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycero-3-PE	5
28154	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycero-3-PC	10
27591	1-Palmitoyl-d ₉ -2-Palmitoyl- <i>sn</i> -glycerol	5
27592	1-Palmitoyl-d ₉ -2,3-Dipalmitoyl- <i>sn</i> -glycerol	10
28123	Cholesteryl Palmitate-d ₉	100

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