

PRODUCT INFORMATION



C26:1 Sphingomyelin-d₅ (d18:1/26:1(17Z)-d₅)

Item No. 9003463

Formal Name: (2S,3R,E)-2-((Z)-hexacos-17-enamido-25,25,26,26,26-d₅)-3-hydroxyoctadec-4-en-1-yl (2-(trimethylammonio)ethyl) phosphate

Synonyms: 17Z-Hexacosenoyl Sphingomyelin-d₅, N-17Z-hexacosenoyl-D-erythro-Sphingosylphosphorylcholine-d₅, SM(d18:1/26:1(17Z)-d₅), Sphingomyelin-d₅ (d18:1/26:1(17Z)-d₅)

MF: C₄₉H₉₂D₅N₂O₆P

FW: 846.3

Chemical Purity: ≥98% (C26:1 Sphingomyelin (d18:1/26:1(17Z)))

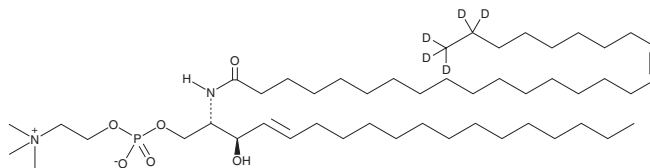
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years



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

Laboratory Procedures

C26:1 Sphingomyelin-d₅ is intended for use as an internal standard for the quantification of C26:1 sphingomyelin by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Description

C26:1 Sphingomyelin is a naturally occurring sphingolipid that has been found in human aqueous humor.¹

Reference

1. Aljohani, A.J., Munguba, G.C., Guerra, Y., et al. Sphingolipids and ceramides in human aqueous humor. *Mol. Vis.* **19**, 1966-1984 (2013).

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, 12/21/2022

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