

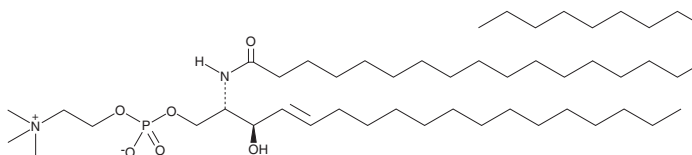
PRODUCT INFORMATION



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

Item No. 9003459

CAS Registry No.: 1448012-56-3
Formal Name: (7S,25Z)-4-hydroxy-7-[(1R,2E)-1-hydroxy-2-hexadecen-1-yl]-N,N,N-trimethyl-9-oxo-3,5-dioxo-8-aza-4-phosphatetriacont-25-en-1-aminium, 4-oxide, inner salt
Synonyms: 17Z-Hexacosenoyl Sphingomyelin, N-(17Z-hexacosenoyl)-D-erythro-Sphingosylphosphorylcholine, SM(d18:1/26:1(17Z)), Sphingomyelin (d18:1/26:1(17Z))



MF: C₄₉H₉₇N₂O₆P
FW: 841.3
Purity: ≥98%
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.

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, 11/02/2022

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