

PRODUCT INFORMATION



C2 Sphingomyelin (d18:1/2:0)

Item No. 24353

CAS Registry No.: 74713-54-5
Formal Name: 4-hydroxy-7-(1-hydroxy-2-hexadecen-1-yl)-N,N,N-trimethyl-9-oxo-3,5-dioxo-8-aza-4-phosphadecan-1-aminium, 4-oxide, inner salt

Synonyms: Acetyl Sphingomyelin, N-Acetyl Sphingosylphosphorylcholine, SM(d18:1/2:0), Sphingomyelin (d18:1/2:0)

MF: C₂₅H₅₁N₂O₆P

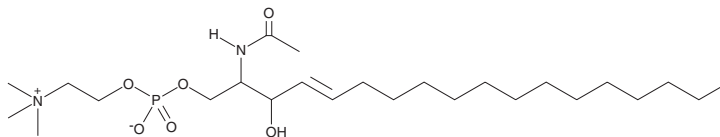
FW: 506.7

Purity: ≥98%

Supplied as: A 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

C2 Sphingomyelin (d18:1/2:0) is supplied as a solid. A stock solution may be made by dissolving the C2 sphingomyelin (d18:1/2:0) in the solvent of choice, which should be purged with an inert gas. C2 Sphingomyelin (d18:1/2:0) is soluble in ethanol and a 2:1 solution of chloroform:methanol.

Description

C2 Sphingomyelin is a synthetic cell-permeable derivative of sphingomyelin (Item Nos. 22674 | 24345). It is composed of a sphingomyelin backbone amine-linked to a C2 fatty acid chain. This product is a mixture of the D-erythro and L-threo isomers of C2 sphingomyelin. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

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, 11/07/2022

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