

# PRODUCT INFORMATION



## C26 Sphingomyelin (d18:1/26:0)

Item No. 9003457

CAS Registry No.: 221097-57-0

Formal Name: (7S)-4-hydroxy-7-[(1R,2E)-1-hydroxy-2-hexadecen-1-yl]-N,N,N-trimethyl-9-oxo-3,5-dioxo-8-aza-4-phosphatetriacontan-1-aminium, 4-oxide, inner salt

Synonyms: Ceroyl Sphingomyelin, N-Hexacosanoyl-D-erythro-Sphingophosphorylcholine, SM(d18:1/26:0), Sphingomyelin (d18:1/26:0)

MF: C<sub>49</sub>H<sub>99</sub>N<sub>2</sub>O<sub>6</sub>P

FW: 843.3

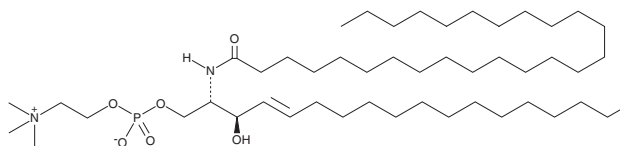
Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years

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



### Description

C26 Sphingomyelin is a sphingolipid that has been found in serum and cell or brain extracts.<sup>1,2,3</sup> C26 sphingomyelin levels are increased in the serum of patients with hemophagocytic lymphohistiocytosis, an inflammatory condition characterized by excessive immune activation.<sup>3</sup> C26 sphingomyelin has been used a component of monolayers to study the influence of sphingomyelin acyl chain length on sphingomyelin-sterol interactions.<sup>4</sup>

### References

1. Wang, J.-R., Zhang, H., Yau, L.D., *et al.* Improved sphingolipidomic approach based on ultra-high performance liquid chromatography and multiple mass spectrometries with application to cellular neurotoxicity. *Anal. Chem.* **86(12)**, 5688-5696 (2014).
2. Willmann, J., Mahlstedt, K., Leibfritz, D., *et al.* Characterization of sphingomyelins in lipid extracts using a HPLC-MS-offline-NMR method. *Anal. Chem.* **79(11)**, 4188-4191 (2007).
3. Jenkins, R.W., Clarke, C.J., Lucas, J.T., Jr., *et al.* Evaluation of the role of secretory sphingomyelinase and bioactive sphingolipids as biomarkers in hemophagocytic lymphohistiocytosis. *Am. J. Hematol.* **88(11)**, E265-E272 (2013).
4. Li, X.M., Momsen, M.M., Brockman, H.L., *et al.* Sterol structure and sphingomyelin acyl chain length modulate lateral packing elasticity and detergent solubility in model membranes. *Biophys. J.* **85(6)**, 3788-801 (2003).

#### 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, 05/05/2020

#### 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