

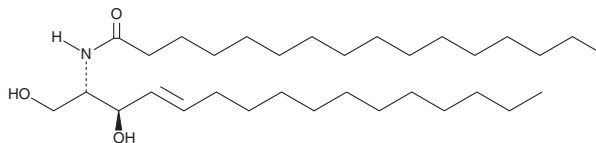
PRODUCT INFORMATION



C16 Ceramide (d16:1/16:0)

Item No. 24426

CAS Registry No.: 123065-37-2
Formal Name: N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-pentadecen-1-yl]-hexadecanamide
Synonyms: Cer(d16:1/16:0), Ceramide (d16:1/16:0), N-hexadecanoyl-D-erythro-Sphingosine C16
MF: $C_{32}H_{63}NO_3$
FW: 509.9
Purity: $\geq 98\%$
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

C16 Ceramide (d16:1/16:0) is supplied as a solid. A stock solution may be made by dissolving the C16 ceramide (d16:1/16:0) in the solvent of choice. C16 Ceramide (d16:1/16:0) is soluble in ethanol (warm), methanol (warm), and chloroform.

Description

C16 Ceramide (d16:1/16:0) is a bioactive sphingolipid that induces ordered- and gel-phase formation and thermal stabilization of palmitoylsphingomyelin-rich domains in palmitoylphosphocholine (POPC) bilayers.¹ It has been found as a component of sphingomyelins (Item Nos. 22674 | 24345) in human erythrocytes.²

References

1. Maule, T., Artetxe, I., Grandell, P.M., *et al.* Importance of the sphingoid base length for the membrane properties of ceramides. *Biophys. J.* **103**(9), 1870-1879 (2012).
2. Myher, J.J., Kuksis, A., and Pind, S. Molecular species of glycerophospholipids and sphingomyelins of human erythrocytes: Improved method of analysis. *Lipids* **24**(5), 396-407 (1989).

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, 12/12/2022

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