

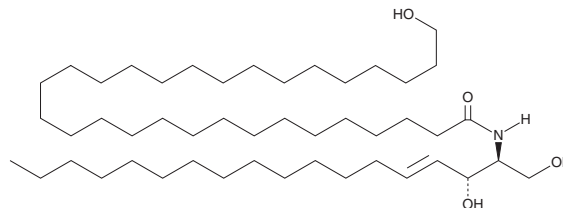
PRODUCT INFORMATION



C30 (ω -hydroxy) Ceramide (d18:1/30:0)

Item No. 22688

CAS Registry No.: 457100-08-2
Formal Name: 30-hydroxy-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-triacontanamide
Synonyms: (ω -hydroxy) Ceramide (d18:1/30:0), (ω -hydroxy) Cer(d18:1/30:0), ω -hydroxy-C30 Sphingosine (d18:1/30:0), ω -OH C30 Ceramide (d18:1/30:0)
MF: C₄₈H₉₅NO₄
FW: 750.3
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

C30 (ω -hydroxy) Ceramide (d18:1/30:0) is supplied as a solid. A stock solution may be made by dissolving the C30 (ω -hydroxy) ceramide (d18:1/30:0) in the solvent of choice. C30 (ω -hydroxy) Ceramide (d18:1/30:0) is soluble in a 5:1 solution of chloroform:methanol. C30 (ω -hydroxy) Ceramide (d18:1/30:0) is also soluble in dimethyl formamide and ethanol (heated).

Description

C30 (ω -hydroxy) Ceramide is a bioactive sphingolipid found in the stratum corneum layer of mammalian epidermis.¹ C30 (ω -hydroxy) Ceramide is vital to skin barrier functions and has a role in maturation of epidermis cells of the spermatozoa/testes. C30 (ω -hydroxy) Ceramide deficiency causes defective skin-water permeability barrier function and neonatal lethality in elongation of very long chain fatty acids-4 (ELOVL4) deficient mice.²

References

1. Sandhoff, R. Very long chain sphingolipids: Tissue expression, function and synthesis. *FEBS Lett.* **584(9)**, 1907-1913 (2010).
2. Li, W., Sandhoff, R., Kono, M., *et al.* Depletion of ceramides with very long chain fatty acids causes defective skin permeability barrier function, and neonatal lethality in ELOVL4 deficient mice. *Int. J. Biol. Sci.* **3(2)**, 120-128 (2007).

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, 01/28/2025

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