

PRODUCT INFORMATION



C32 Ceramide (d18:1/32:0)

Item No. 24424

CAS Registry No.: 34227-73-1
Formal Name: N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-dotriacontanamide

Synonyms: N-dotriacontanoyl-D-erythro-Sphingosine, Cer(d18:1/32:0), Ceramide (d18:1/32:0)

MF: C₅₀H₉₉NO₃

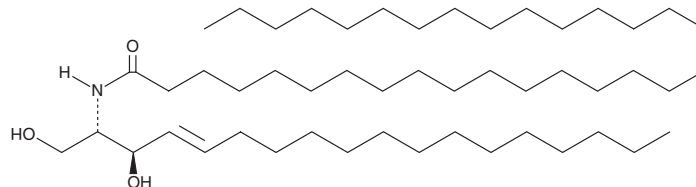
FW: 762.3

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

C32 Ceramide (d18:1/32:0) is supplied as a solid. A stock solution may be made by dissolving the C32 ceramide (d18:1/32:0) in the solvent of choice. C32 Ceramide (d18:1/32:0) is soluble in methanol (warmed), ethanol, (warmed), DMSO (warmed), and in a 2:1 solution of chloroform:methanol.

Description

C32 Ceramide is a naturally occurring ceramide that has been found in the stratum corneum of human and nude mouse skin.^{1,2}

References

1. Farwanah, H., Wohlrab, J., Neubert, R.H., *et al.* Profiling of human stratum corneum ceramides by means of normal phase LC/APCI-MS. *Anal. Bioanal. Chem.* **383**(4), 632-637 (2005).
2. Liou, Y.B., Sheu, M.T., Liu, D.Z., *et al.* Quantitation of ceramides in nude mouse skin by normal-phase liquid chromatography and atmospheric pressure chemical ionization mass spectrometry. *Anal. Biochem.* **401**(1), 107-113 (2010).

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, 12/19/2022

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