

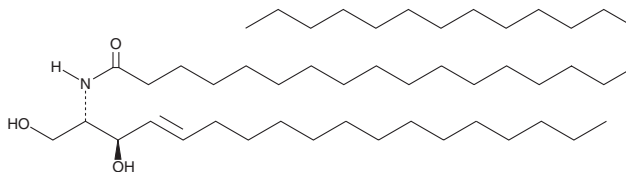
# PRODUCT INFORMATION



## C30 Ceramide (d18:1/30:0)

Item No. 24400

**CAS Registry No.:** 871540-97-5  
**Formal Name:** N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-triacontanamide  
**Synonyms:** Cer(d18:1/30:0), Ceramide (d18:1/30:0), N-triacontanoyl-D-erythro-Sphingosine  
**MF:** C<sub>48</sub>H<sub>95</sub>NO<sub>3</sub>  
**FW:** 734.3  
**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

C30 Ceramide (d18:1/30:0) is supplied as a solid. A stock solution may be made by dissolving the C30 ceramide (d18:1/30:0) in the solvent of choice, which should be purged with an inert gas. C30 Ceramide (d18:1/30:0) is soluble in chloroform and ethanol, and a 2:1 solution of chloroform:methanol.

### Description

C30 Ceramide is a naturally occurring ceramide that has been found in the stratum corneum of human and nude mouse skin.<sup>1,2</sup> C30 Ceramide levels are decreased in the skin of patients with lesional and non-lesional atopic dermatitis.<sup>3</sup>

### 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).
3. Joo, J.-M., Hwang, J.-H., Bae, S., *et al.* Relationship of ceramide-, and free fatty acid-cholesterol ratios in the stratum corneum with skin barrier function of normal, atopic dermatitis lesional and non-lesional skins. *J. Dermatol. Sci.* **77**(1), 71-74 (2015).

#### 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, 11/16/2022

#### 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