

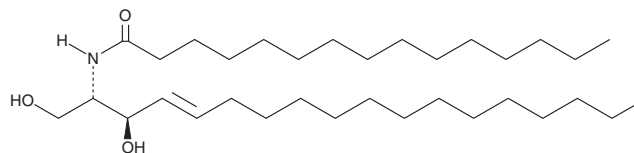
# PRODUCT INFORMATION



## C15 Ceramide (d18:1/15:0)

Item No. 24393

**CAS Registry No.:** 67492-15-3  
**Formal Name:** N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-pentadecanamide  
**Synonyms:** Cer(d18:1/15:0), Ceramide (d18:1/15:0), N-pentadecanoyl-D-erythro-Sphingosine  
**MF:** C<sub>33</sub>H<sub>65</sub>NO<sub>3</sub>  
**FW:** 523.9  
**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

C15 Ceramide (d18:1/15:0) is supplied as a solid. A stock solution may be made by dissolving the C15 ceramide (d18:1/15:0) in the solvent of choice. C15 Ceramide (d18:1/15:0) is soluble in organic solvents such as chloroform, ethanol (warmed), and methanol (warmed).

### Description

C15 Ceramide is a naturally occurring ceramide that has been found in *L. nidifica* and *A. glomerata* extracts, human erythrocytes, and the stratum corneum of *P. domesticus*.<sup>1-3</sup>

### References

- Muñoz-García, A., Ro, J., Brown, J.C., *et al.* Identification of complex mixtures of sphingolipids in the stratum corneum by reversed-phase high-performance liquid chromatography and atmospheric pressure photospray ionization mass spectrometry. *J. Chromatogr. A.* **1133(1-2)**, 58-68 (2006).
- 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).
- Cardellina, J.H. and Moore, R.E. Sphingosine derivatives from red algae of the ceramiales. *Phytochemistry* **17(3)**, 554-555 (1978).

#### 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