

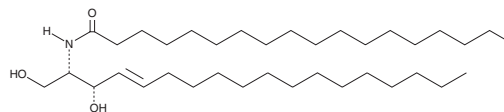
# PRODUCT INFORMATION



## C18 L-threo Ceramide (d18:1/18:0)

Item No. 24399

**CAS Registry No.:** 95037-06-2  
**Formal Name:** N-[(1S,2S,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-octadecanamide  
**Synonyms:** N-octadecanoyl-L-threo-Sphingosine, L-threo Cer(d18:1/18:0), L-threo Ceramide (d18:1/18:0)  
**MF:** C<sub>36</sub>H<sub>71</sub>NO<sub>3</sub>  
**FW:** 566.0  
**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

C18 L-threo Ceramide (d18:1/18:0) is supplied as a solid. A stock solution may be made by dissolving the C18 L-threo ceramide (d18:1/18:0) in the solvent of choice, which should be purged with an inert gas. C18 L-threo Ceramide (d18:1/18:0) is soluble in a 90:10 solution of chloroform:methanol.

### Description

C18 L-threo Ceramide is a synthetic ceramide and stereoisomer of C18 ceramide (Item No. 19556) that has been used for structural characterization of natural sphingolipids.<sup>1</sup> It inhibits rat brain mitochondrial ceramidase (mt-CDase) with an IC<sub>50</sub> value of 8.8 μM.<sup>2</sup>

### References

1. Sarmientos, F., Schwarzmann, G., and Sandhoff, K. Direct evidence by carbon-13 NMR spectroscopy for the erythro configuration of the sphingoid moiety in Gaucher cerebroside and other natural sphingolipids. *Eur. J. Biochem.* **146**(1), 59-64 (1985).
2. Usta, J., El Bawab, S., Roddy, P., et al. Structural requirements of ceramide and sphingosine based inhibitors of mitochondrial ceramidase. *Biochemistry* **40**(32), 9657-9668 (2001).

#### 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, 12/17/2022

#### 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](http://WWW.CAYMANCHEM.COM)