

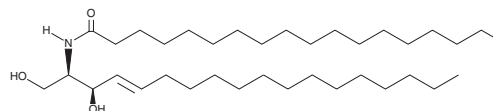
PRODUCT INFORMATION



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

Item No. 33676

CAS Registry No.: 252039-53-5
Formal Name: N-[(1R,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-octadecanamide
Synonyms: D-threo Cer(d18:1/18:0),
D-threo Ceramide (d18:1/18:0),
N-octadecanoyl-D-threo-Sphingosine,
N-stearoyl-D-threo-Sphingosine
MF: C₃₆H₇₁NO₃
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 D-threo Ceramide (d18:1/18:0) is supplied as a solid. A stock solution may be made by dissolving the C18 D-threo ceramide (d18:1/18:0) in the solvent of choice, which should be purged with an inert gas. C18 D-threo Ceramide (d18:1/18:0) is soluble in ethanol, chloroform, DMSO, and dimethyl formamide.

Description

C18 D-threo Ceramide (d18:1/18:0) is a synthetic ceramide and stereoisomer of C18 ceramide (Item No. 19556).¹ It inhibits rat brain mitochondrial ceramidase (mt-CDase) with an IC₅₀ value of 16.6 μM.²

References

1. Koolath, S., Murai, Y., Suga, Y., *et al.* Chiral combinatorial preparation and biological evaluation of unique ceramides for inhibition of sphingomyelin synthase. *Chirality* **32(3)**, 308-313 (2019).
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, 11/15/2022

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