

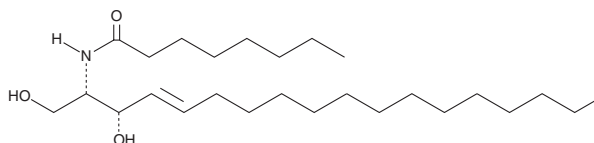
PRODUCT INFORMATION



C8 L-threo Ceramide (d18:1/8:0)

Item No. 24392

CAS Registry No.: 175892-44-1
Formal Name: N-[(1S,2S,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-octanamide
Synonyms: L-threo Cer(d18:1/8:0), L-threo Ceramide (d18:1/8:0), N-octanoyl-L-threo-Sphingosine
MF: C₂₆H₅₁NO₃
FW: 425.7
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

C8 L-threo Ceramide (d18:1/8:0) is supplied as a solid. A stock solution may be made by dissolving the C8 L-threo ceramide (d18:1/8:0) in the solvent of choice, which should be purged with an inert gas. C8 L-threo ceramide (d18:1/8:0) is soluble in chloroform, ethanol, DMSO, and dimethyl formamide (DMF). The solubility of C8 L-threo ceramide (d18:1/8:0) in DMF is approximately C8 L-threo ceramide 5 mg/ml.

Description

C8 L-threo Ceramide is a bioactive sphingolipid and cell-permeable analog of naturally occurring ceramides.¹ It is cytotoxic to U937 cells (IC₅₀ = 10 μM) and, like C8 D-threo ceramide (Item No. 24391), induces nuclear DNA fragmentation 5- to 6-fold more potently than C8 D-erythro ceramide (Item No. 62540).^{1,2} C8 L-threo Ceramide also enhances *V. cholerae* cytolysin pore formation in liposome lipid membranes, as measured by calcein release, with a 50% release dose (RD₅₀) value of ~30 μg/ml.³

References

1. Karasavvas, N., Erukulla, R.K., Bittman, R., *et al.* Stereospecific induction of apoptosis in U937 cells by N-octanoyl-sphingosine stereoisomers and N-octyl-sphingosine. The ceramide amide group is not required for apoptosis. *Eur. J. Biochem.* **236**(2), 729-737 (1996).
2. Chang, Y.-T., Choi, J., Ding, S., *et al.* The synthesis and biological characterization of a ceramide library. *J. Am. Chem. Soc.* **124**(9), 1856-1857 (2002).
3. Zitzer, A., Bittman, R., Verbicky, C.A., *et al.* Coupling of cholesterol and cone-shaped lipids in bilayers augments membrane permeabilization by the cholesterol-specific toxins streptolysin O and *Vibrio cholerae* cytolysin. *J. Biol. Chem.* **276**(18), 14628-14633 (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, 06/26/2024

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