

PRODUCT INFORMATION



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

Item No. 24398

CAS Registry No.: 252039-52-4
Formal Name: N-[(1R,2S,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-octadecanamide

Synonyms: L-erythro Cer(d18:1/18:0), L-erythro Ceramide (d18:1/18:0), N-octadecanoyl-L-erythro-Sphingosine

MF: C₃₆H₇₁NO₃

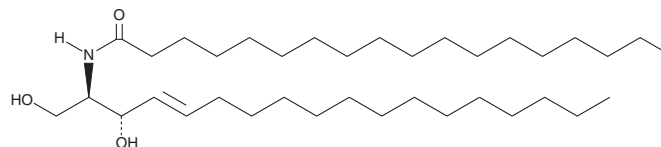
FW: 566.0

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 L-erythro Ceramide (d18:1/18:0) is supplied as a solid. A stock solution may be made by dissolving the C18 L-erythro ceramide (d18:1/18:0) in the solvent of choice, which should be purged with an inert gas. C18 L-erythro Ceramide (d18:1/18:0) is soluble in organic solvents such as chloroform, ethanol, and DMSO.

Description

C18 L-erythro Ceramide is a naturally occurring ceramide and stereoisomer of C18 ceramide (Item No. 19556).¹ It inhibits rat brain mitochondrial ceramidase (mt-CDase) with an IC₅₀ value of 0.11 mol% (8.8 μM).

Reference

1. 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, 04/09/2024

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