

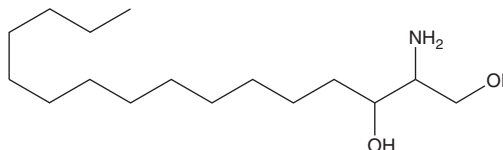
PRODUCT INFORMATION



DL-erythro/threo Sphinganine (d16:0)

Item No. 24376

CAS Registry No.: 50731-08-3
Formal Name: 2-amino-1,3-hexadecanediol
Synonym: DL-erythro/threo Sphinganine C16
MF: $C_{16}H_{35}NO_2$
FW: 273.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

DL-erythro/threo Sphinganine (d16:0) is supplied as a solid. A stock solution may be made by dissolving the DL-erythro/threo sphinganine (d16:0) in the solvent of choice. DL-erythro/threo Sphinganine (d16:0) is soluble in organic solvents such as chloroform, methanol, ethanol, and DMSO, which should be purged with an inert gas.

Description

DL-erythro/threo Sphinganine (d16:0) is a sphingolipid that is decreased in rats following long-term, low-dose administration of dimethoate and is used as a biomarker for dimethoate exposure.¹ It has been found in the fermentation broth of *Bacterium SRCnm*, which has antimicrobial and antioxidant activities.² This product is a mixture of sphinganine (d16:0), L-erythro sphinganine (d16:0), D-threo sphinganine (d16:0), and L-threo sphinganine (d16:0).

References

1. Feng, Z., Sun, X., Yang, J., et al. Metabonomics analysis of urine and plasma from rats given long-term and low-dose dimethoate by ultra-performance liquid chromatography-mass spectrometry. *Chem. Biol. Interact.* **199(3)**, 143-153 (2012).
2. Ahmed, E.F., Hassan, H.M., Rateb, M.E., et al. A comparative biochemical study on two marine endophytes, *Bacterium SRCnm* and *Bacillus sp. JS*, isolated from red sea algae. *Pak. J. Pharm. Sci.* **29(1)**, 17-26 (2016).

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