

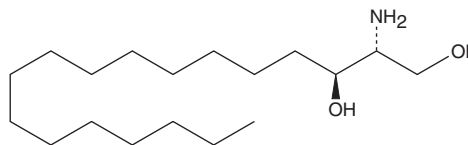
PRODUCT INFORMATION



L-erythro Sphinganine (d18:0)

Item No. 24374

CAS Registry No.: 6036-76-6
Formal Name: 2R-amino-1,3S-octadecanediol
Synonyms: L-erythro Dihydrosphingosine,
C18 L-erythro Sphinganine (d18:0)
MF: C₁₈H₃₉NO₂
FW: 301.5
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

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

Description

L-erythro Sphinganine (d18:0) is a synthetic bioactive sphingolipid and stereoisomer of sphinganine (d18:0) (Item No. 10007945) and D-threo sphinganine (d18:0) (Item No. 24375).¹ It induces autophagy in HCT116 cells when used at a concentration of 12 μM. L-erythro Sphinganine induces cell cycle arrest of 7R4-LCB1 yeast cells.¹ It is metabolized via sphinganine N-acyltransferase and sphinganine kinase *in vivo* in rat liver.²

References

1. Coward, J., Ambrosini, G., Musi, E., *et al.* Safingol (L-threo-sphinganine) induces autophagy in solid tumor cells through inhibition of PKC and the PI3-kinase pathway. *Autophagy* **5(2)**, 184-193 (2009).
2. Stoffel, W. and Bister, K. Stereospecificities in the metabolic reactions of the four isomeric sphinganine (dihydrosphingosines) in rat liver. *Hoppe-Seylers Z. Physiol. Chem.* **354(2)**, 169-181 (1973).

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, 10/04/2022

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