

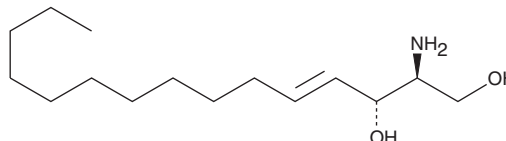
PRODUCT INFORMATION



Sphingosine (d15:1)

Item No. 10007901

CAS Registry No.: 86555-28-4
Formal Name: 2S-amino-4E-pentadecene-1,3R-diol
Synonym: D-erythro-Sphingosine C-15
MF: $C_{15}H_{31}NO_2$
FW: 257.4
Purity: $\geq 98\%$
Supplied as: A crystalline 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

Sphingosine (d15:1) is supplied as a crystalline solid. A stock solution may be made by dissolving the sphingosine (d15:1) in the solvent of choice, which should be purged with an inert gas. Sphingosine (d15:1) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of sphingosine (d15:1) is miscible in ethanol, whereas the solubility is approximately 2 and 10 mg/ml in DMSO and DMF, respectively.

Description

Sphingosine is an amino alcohol most commonly characterized by an 18-carbon unsaturated hydrocarbon chain (Sphingosine (d18:1), Item No. 10007907). However, the hydrocarbon chain length of sphingosine, and the related dihydrosphingosine, can vary from 12-26 carbons in mammalian tissues.^{1,2} Sphingosine (d15:1) is a naturally-occurring but rare form of sphingosine. It is intended to be used as an internal standard in the analysis of sphingoid compounds by chromatographic or spectrometric methods.

References

1. Stewart, M.E. and Downing, D.T. Free sphingosines of human skin include 6-hydroxysphingosine and unusually long-chain dihydrosphingosines. *J. Invest. Dermatol.* **105**, 613-618 (1995).
2. Pruett, S.T., Bushnev, A., Hagedorn, K., et al. Biodiversity of sphingoid bases ("sphingosines") and related amino alcohols. *J. Lipid Res.* **49**, 1621-1639 (2008).

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, 10/05/2022

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