

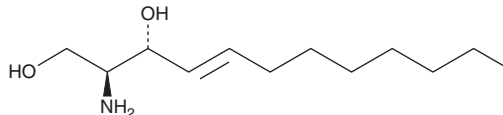
PRODUCT INFORMATION



Sphingosine (d12:1)

Item No. 26893

CAS Registry No.: 128427-86-1
Formal Name: 2S-amino-4E-dodecene-1,3R-diol
Synonym: D-erythro-Sphingosine C12
MF: C₁₂H₂₅NO₂
FW: 215.3
Purity: ≥95%
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 (d12:1) is supplied as a crystalline solid. A stock solution may be made by dissolving the sphingosine (d12:1) in the solvent of choice. Sphingosine (d12:1) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of sphingosine (d12:1) in these solvents is approximately 2 and 10 mg/ml, respectively. Sphingosine (d12:1) is also miscible in ethanol.

Description

Sphingosine (d12:1) is a short-chain sphingolipid.^{1,2} It decreases serine palmitoyltransferase activity in primary cultured mouse cerebellar cells in a concentration-dependent manner.¹ Sphingosine (d12:1) (0.5-50 μM) decreases production of sphingosine (d18:1) (Item No. 10007907), lactosylceramide (Item Nos. 27197 | 16983), galactosylceramide (Item No. 24322), and a variety of gangliosides in mouse cerebellar granule cells.²

References

1. Mandon, E.C., van Echten, G., Birk, R., *et al.* Sphingolipid biosynthesis in cultured neurons: Down-regulation of serine palmitoyltransferase by sphingoid bases. *Eur. J. Biochem.* **198(3)**, 667-674 (1991).
2. van Echten, G., Birk, R., Brenner-Weiss, G., *et al.* Modulation of sphingolipid biosynthesis in primary cultured neurons by long chain bases. *J. Biol. Chem.* **265(16)**, 9333-9339 (1990).

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, 12/13/2022

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