

PRODUCT INFORMATION



Sphingosine (d18:1(14Z))

Item No. 27019

Formal Name: (2S,3R,Z)-2-amino-octadec-14-ene-1,3-diol

Synonym: Sphing-14Z-ene

MF: C₁₈H₃₇NO₂

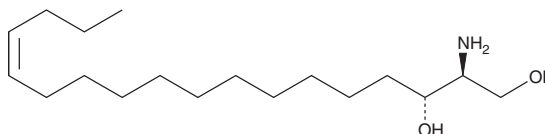
FW: 299.5

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 (d18:1(14Z)) is supplied as a crystalline solid. A stock solution may be made by dissolving the sphingosine (d18:1(14Z)) in the solvent of choice, which should be purged with an inert gas. Sphingosine (d18:1(14Z)) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of sphingosine (d18:1(14Z)) in these solvents is approximately 2 and 10 mg/ml, respectively. Sphingosine (d18:1(14Z)) is also miscible in ethanol.

Description

Sphingosine (d18:1(14Z)) is an atypical sphingolipid that contains a *cis* double bond at the 14-15 position rather than a *trans* double bond at the 4-5 position as in sphingosine (d18:1) (Item No. 10007907).

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, 02/16/2024

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