

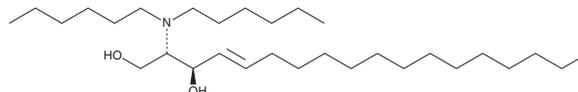
PRODUCT INFORMATION



N,N-dihexyl Sphingosine (d18:1/6:0/6:0)

Item No. 24385

CAS Registry No.: 2377379-53-6
Formal name: (2S,3R,E)-2-(dihexylamino)octadec-4-ene-1,3-diol
Synonym: N,N-dihexyl-D-erythro-Sphingosine
MF: C₃₀H₆₁NO₂
FW: 467.8
Purity: ≥95%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

N,N-dihexyl Sphingosine (d18:1/6:0/6:0) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. N,N-dihexyl Sphingosine (d18:1/6:0/6:0) is soluble in organic solvents such as chloroform, and methanol, which should be purged with an inert gas.

Description

N,N-dihexyl Sphingosine is a synthetic analog of C18 ceramide (Item No. 19556) that has two hexyl groups acylated to sphingosine in place of the naturally occurring stearoyl fatty acid.

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, 01/06/2023

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