

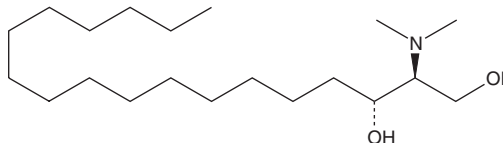
PRODUCT INFORMATION



N,N-dimethyl Sphinganine (d18:0)

Item No. 22504

CAS Registry No.: 17267-46-8
Formal Name: 2S-(dimethylamino)-1,3R-octadecanediol
Synonym: N,N-dimethyl-D-erythro-Dihydrosphingosine
MF: C₂₀H₄₃NO₂
FW: 329.6
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-dimethyl Sphinganine (d18:0) is supplied as a solution in ethanol. To change the solvent, simply evaporate the N,N-dimethyl sphinganine (d18:0) under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of N,N-dimethyl sphinganine (d18:0) in these solvents is approximately 2 and 10 mg/ml, respectively.

Description

N,N-dimethyl Sphinganine (d18:0) is a sphingolipid that increases during development of *L. donovani* promastigotes.¹

Reference

1. Silva, A.M., Cordeiro-da-Silva, A., and Coombs, G.H. Metabolic variation during development in culture of *Leishmania donovani* promastigotes. *PLoS Negl. Trop. Dis.* **5(12)**, e1451 (2011).

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, 04/20/2018

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