

PRODUCT INFORMATION



C6 NBD L-threo dihydro Ceramide (d18:0/6:0)

Item No. 24440

Formal Name: N-(1,3R-dihydroxyoctadecan-2R-yl)-6-((7-nitrobenzo[c][1,2,5]oxadiazol-4-yl)amino)hexanamide

Synonyms: N-hexanoyl-NBD L-threo-Dihydrosphingosine, NBD L-threo Cer(d18:0/6:0), NBD L-threo Ceramide (d18:0/6:0)

MF: C₃₀H₅₁N₅O₆

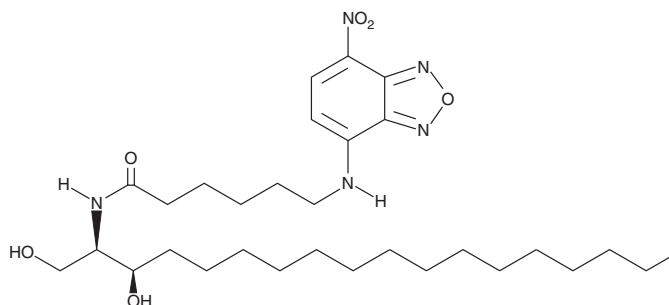
FW: 577.8

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

C6 NBD L-threo dihydro ceramide (d18:0/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 NBD L-threo dihydro ceramide (d18:0/6:0) in the solvent of choice. C6 NBD L-threo dihydro ceramide (d18:0/6:0) is soluble in methanol and in a 2:1 solution of chloroform:methanol, which should be purged with an inert gas..

Description

C6 NBD L-threo dihydro ceramide is a sphingolipid that has been tagged with a fluorescent C6 nitrobenzoxadiazole (C6 NBD; Item No. 18871) group. It is an isomer of C6 NBD dihydro ceramide (Item No. 10007957).

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, 02/09/2018

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