

PRODUCT INFORMATION

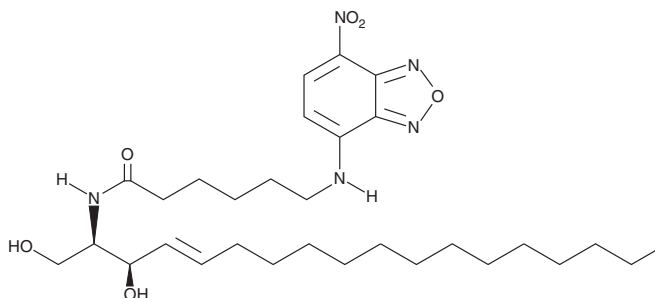


C6 NBD L-threo Ceramide (d18:1/6:0)

Item No. 24439

CAS Registry No.: 114301-95-0
Formal Name: [R-[R*,R*-(E)]]-N-[2-hydroxy-1-(hydroxymethyl)-3-heptadecenyl]-6-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-hexanamide
Synonyms: NBD L-threo Cer(d18:1/6:0), NBD L-threo Ceramide (d18:1/6:0), N-hexanoyl-NBD L-threo-Sphingosine

MF: C₃₀H₄₉N₅O₆
FW: 575.7
Purity: ≥98%
Supplied as: A 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

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

Description

C6 NBD L-threo ceramide is a biologically active and cell-permeable analog of naturally occurring ceramides that is tagged with a fluorescent C6 nitrobenzoxadiazole (C6 NBD; Item No. 18871) group. It is rapidly transferred between liposomes, labels the Golgi apparatus, and is metabolized to C6 NBD sphingomyelin (Item No. 24328) in BHK cells and V79 fibroblasts.¹

Reference

1. Pagano, R.E. and Martin, O.C. A series of fluorescent N-acylsphingosines: Synthesis, physical properties, and studies in cultured cells. *Biochemistry* **27**(12), 4439-4445 (1988).

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, 12/17/2022

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