

PRODUCT INFORMATION



C6 NBD Phytoceramide (t18:0/6:0)

Item No. 24326

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

Synonyms: C6 NBD Phytosphingosine (t18:0/6:0), N-hexanoyl-NBD Phytosphingosine, NBD Cer(t18:0/6:0), NBD Ceramide (t18:0/6:0)

MF: C₃₀H₅₁N₅O₇

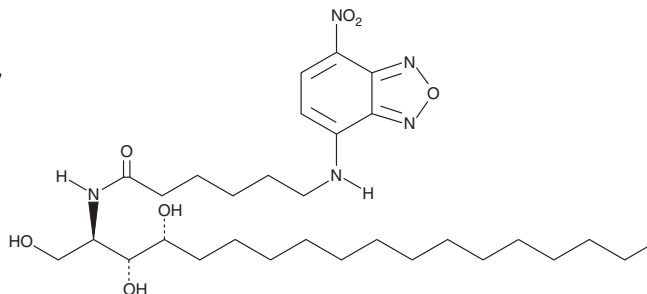
FW: 593.8

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 phytoceramide (t18:0/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 NBD phytoceramide (t18:0/6:0) in the solvent of choice, which should be purged with an inert gas. C6 NBD phytoceramide (t18:0/6:0) is soluble in the organic solvent methanol and chloroform/methanol (2:1).

Description

C6 NBD phytoceramide is a biologically active derivative of C6 phytoceramide (Item No. 24438) that is tagged with a fluorescent C6 nitrobenzoxadiazole (C6 NBD; Item No. 18871) group.¹ C6 NBD phytoceramide localizes to the Golgi apparatus of cultured fibroblasts and is metabolized to fluorescent sphingomyelin and glucosylceramide.

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, 11/16/2022

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