

PRODUCT INFORMATION



C12 NBD Phytoceramide (t18:0/12:0)

Item No. 24327

Formal Name: 12-((7-nitrobenzo[c][1,2,5]oxadiazol-4-yl)amino)-N-((2R,3R,4R)-1,3,4-trihydroxyoctadecan-2-yl)dodecanamide

Synonyms: C12 NBD Phytosphingosine (t18:0/12:0), NBD Cer(t18:0/12:0), NBD Ceramide (t18:0/12:0), N-dodecanoyl-NBD Phytosphingosine

MF: C₃₆H₆₃N₅O₇

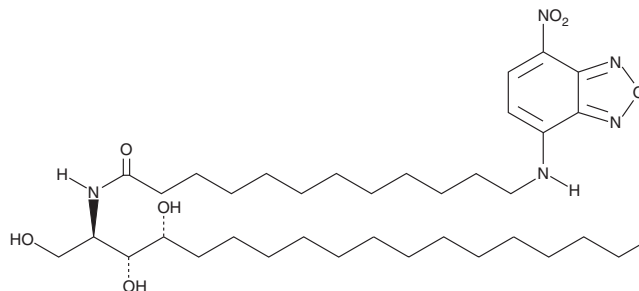
FW: 677.9

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

C12 NBD phytoceramide (t18:0/12:0) is supplied as a solid. A stock solution may be made by dissolving the C12 NBD phytoceramide (t18:0/12:0) in the solvent of choice, which should be purged with an inert gas. C12 NBD phytoceramide (t18:0/12:0) is soluble in methanol and a 2:1 solution of chloroform:methanol.

Description

C12 NBD phytoceramide is a biologically active derivative of phytosphingosine (Item No. 20217) that is tagged with a fluorescent C12 nitrobenzoxadiazole (C12 NBD) group. C12 NBD phytosphingosine has been used to study the metabolism of ceramides using microsomes expressing the *S. cerevisiae* diacylglycerol Lro1p.¹

Reference

1. Voynova, N.S., Vionnet, C., Ejsing, C.S., *et al.* A novel pathway of ceramide metabolism in *Saccharomyces cerevisiae*. *Biochem. J.* **447**(1), 103-114 (2012).

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