

PRODUCT INFORMATION



C12 NBD Galactosylceramide (d18:1/12:0)

Item No. 22851

CAS Registry No.: 474942-98-8
Formal Name: N-[(1S,2R,3E)-1-[(β-D-galactopyranosyloxy)methyl]-2-hydroxy-3-heptadecen-1-yl]-12-[[7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-dodecanamide

Synonyms: N-Dodecanoyl-NBD-Galactosylceramide, C12 NBD Galactosylceramide, NBD Galactosylceramide (d18:1/12:0), C12-NBD GalCer, NBD GalCer(d18:1/12:0)

MF: C₄₂H₇₁N₅O₁₁

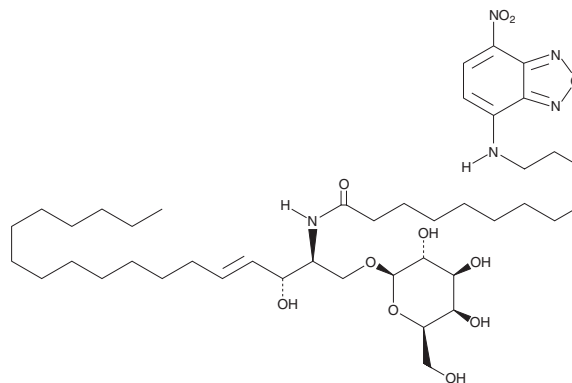
FW: 822.0

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 galactosylceramide (d18:1/12:0) is supplied as a solid. A stock solution may be made by dissolving the C12 NBD galactosylceramide (d18:1/12:0) in the solvent of choice. C12 NBD galactosylceramide (d18:1/12:0) is soluble in chloroform and DMSO and a 2:1 solution of chloroform:methanol.

Description

C12 NBD galactosylceramide is a biologically active derivative of galactosylceramide that is tagged with a fluorescent C12 nitrobenzoxadiazole (C12 NBD) group.¹ C12 NBD galactosylceramide has been used as an internal standard to detect products of C12 NBD ceramide glycosylation. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Reference

1. Turáková, K., Pavlíková, L., Messingerová, L., *et al.* Reduced UDP-glucose levels are associated with P-glycoprotein over-expression in L1210 cells and limit glucosylceramide synthase activity. *Anticancer Res.* **35**(5), 2627-2634 (2015).

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, 11/28/2022

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