

PRODUCT INFORMATION



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

Item No. 22829

CAS Registry No.: 474943-06-1

Formal Name: N-[(1S,2R,3E)-1-[[[4-O-β-D-galactopyranosyl-β-D-glucopyranosyl]oxy]methyl]-2-hydroxy-3-heptadecen-1-yl]-12-[(7-nitro-2,1,3-benzoxadiazol-4-yl)amino]-dodecanamide

Synonyms: C12 NBD Lactosylceramide, NBD LacCer(d18:1/12:0), NBD Lactosylceramide (d18:1/12:0), N-dodecanoyl-NBD-lactosylceramide

MF: C₄₈H₈₁N₅O₁₆

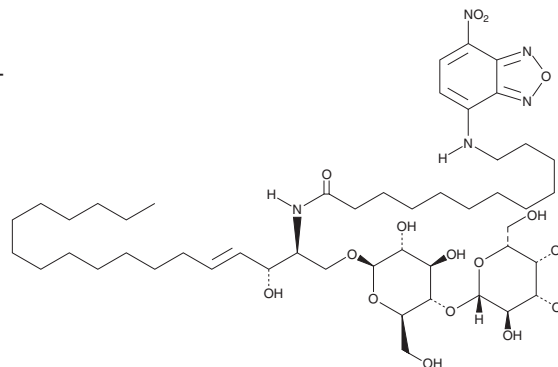
FW: 984.2

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

Description

C12 NBD Lactosylceramide (d18:1/12:0) is a derivative of lactosylceramides (d18:1/acyl mixture) (Item No. 16983) that is tagged with a fluorescent C-12 nitrobenzoxadiazole (C-12 NBD) group. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

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, 12/02/2022

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