

PRODUCT INFORMATION



C12 NBD 3'-sulfo Galactosylceramide (d18:1/12:0)

Item No. 24627

Formal Name: (2R,3R,4S,5S,6R)-3,5-dihydroxy-2-(((2S,3R,E)-3-hydroxy-2-(12-((7-nitrobenzo[c][1,2,5]oxadiazol-4-yl)amino)dodecanamido)octadec-4-en-1-yl)oxy)-6-(hydroxymethyl)tetrahydro-2H-pyran-4-yl hydrogen sulfate

Synonyms: C12 NBD Sulfatide, NBD 3'-sulfo Galactosylceramide (d18:1/12:0), NBD (3'-sulfo)Gal β -Cer(d18:1/12:0), N-dodecanoyl-NBD-Sulfatide

MF: C₄₂H₇₁N₅O₁₄S

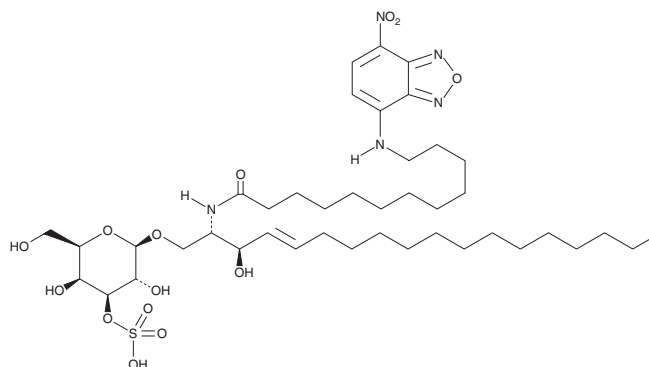
FW: 902.1

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

C12 NBD 3'-sulfo galactosylceramide (d18:1/12:0) is supplied as a solid. A stock solution may be made by dissolving the C12 NBD 3'-sulfo galactosylceramide (d18:1/12:0) in the solvent of choice, which should be purged with an inert gas. C12 NBD 3'-sulfo galactosylceramide (d18:1/12:0) is soluble in a 2:1 solution of chloroform:methanol.

Description

C12 NBD 3'-sulfo galactosylceramide is a sphingolipid that is tagged with a fluorescent C12 nitrobenzoxadiazole (C12 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, 11/10/2022

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