

PRODUCT INFORMATION



C19 3'-sulfo Galactosylceramide (d18:1/19:0)

Item No. 24866

Formal Name: (2R)-3R,5S-dihydroxy-2-(((E)-3R-hydroxy-2S-nonadecanamidoctadec-4-en-1-yl)oxy)-6R-(hydroxymethyl)tetrahydro-2H-pyran-4S-yl hydrogen sulfate

Synonyms: (3'-sulfo)Gal β -Cer(d18:1/19:0), N-Nonadecanoyl Sulfatide, C19 Sulfatide

MF: C₄₃H₈₃NO₁₁S

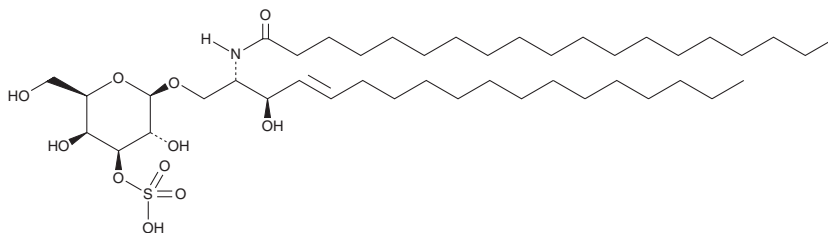
FW: 822.2

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

C19 3'-sulfo Galactosylceramide (d18:1/19:0) is supplied as a solid. A stock solution may be made by dissolving the C19 3'-sulfo galactosylceramide (d18:1/19:0) in the solvent of choice. C19 3'-sulfo Galactosylceramide (d18:1/19:0) is soluble in DMSO, dimethyl formamide, and a 2:1 solution of chloroform:methanol.

Description

C19 3'-sulfo Galactosylceramide is a synthetic derivative of sulfatide (Item No. 24323). 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, 12/19/2022

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