

PRODUCT INFORMATION



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

Item No. 24349

CAS Registry No.: 852100-88-0

Formal Name: N-[(1S,2R,3E)-2-hydroxy-1-[[3-O-sulfo-β-D-galactopyranosyl]oxy]methyl]-3-heptadecen-1-yl]-dodecanamide

Synonyms: N-Dodecanoyl Sulfatide,
(3'-sulfo)Galβ-Cer(d18:1/12:0),
3'-sulfo Galactosylceramide (d18:1/12:0),
C12 Sulfatide

MF: C₃₆H₆₉NO₁₁S

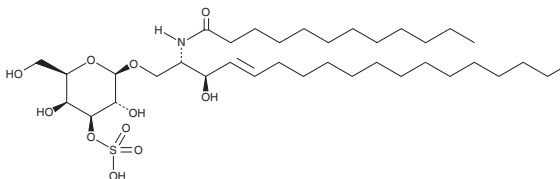
FW: 724.0

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

Description

C12 3'-sulfo Galactosylceramide is a member of the sulfatide class of glycolipids. It has been used as an internal standard for the quantification of sulfatides.¹ As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Reference

1. Wang, J.-R., Zhang, H., Yau, L.D., *et al.* Improved sphingolipidomic approach based on ultra-high performance liquid chromatography and multiple mass spectrometries with application to cellular neurotoxicity. *Anal. Chem.* **86**(12), 5688-5696 (2014).

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, 10/18/2022

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