

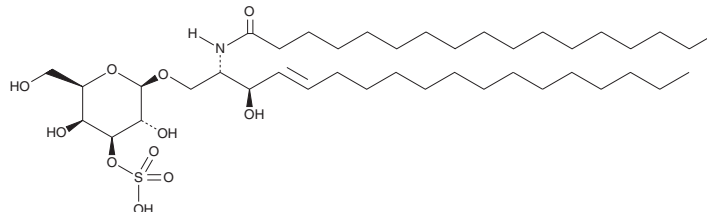
PRODUCT INFORMATION



C17 3'-sulfo Galactosylceramide (d18:1/17:0)

Item No. 24348

CAS Registry No.: 1071654-84-6
Formal Name: N-[(1S,2R,3E)-2-hydroxy-1-[[[3-O-sulfo-β-D-galactopyranosyl]oxy]methyl]-3-heptadecen-1-yl]-heptadecanamide
Synonyms: (3'-sulfo)Galβ-Cer(d18:1/17:0), N-Heptadecanoyl Sulfatide, C17 Sulfatide
MF: C₄₁H₇₉NO₁₁S
FW: 794.1
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

C17 3'-sulfo Galactosylceramide (d18:1/17:0) is supplied as a solid. A stock solution may be made by dissolving the C17 3'-sulfo galactosylceramide (d18:1/17:0) in the solvent of choice. C17 3'-sulfo Galactosylceramide (d18:1/17:0) is soluble in DMSO, dimethyl formamide, and a 2:1 solution of chloroform:methanol.

Description

C17 3'-sulfo Galactosylceramide is a member of the sulfatide class of glycolipids. C17 3'-sulfo Galactosylceramide has been used as an internal standard for the quantification of sulfatides in urine and blood samples from patients with metachromatic leukodystrophy, a disorder characterized by arylsulfatase A deficiency, leading to sulfatide accumulation.^{1,2} As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. Barcenas, M., Suhr, T.R., Scott, C.R., *et al.* Quantification of sulfatides in dried blood and urine spots from metachromatic leukodystrophy patients by liquid chromatography/electrospray tandem mass spectrometry. *Clin. Chim. Acta* **433**, 39-43 (2014).
2. Cui, Y., Colsch, B., Afonso, C., *et al.* Synthetic sulfogalactosylceramide (sulfatide) and its use for the mass spectrometric quantitative urinary determination in metachromatic leukodystrophies. *Glycoconj. J.* **25(2)**, 147-155 (2008).

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, 11/11/2022

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