

PRODUCT INFORMATION



Lyso-globotetraosylceramide (d18:1)

Item No. 29134

Formal Name: N-((2S,3R,4R,5R,6R)-2-(((2R,3R,4S,5S,6R)-2-(((2R,3R,4R,5R,6S)-6-(((2R,3S,4R,5R,6R)-6-(((2S,3R,E)-2-amino-3-hydroxyoctadec-4-en-1-yl)oxy)-4,5-dihydroxy-2-(hydroxymethyl) tetrahydro-2H-pyran-3-yl)oxy)-4,5-dihydroxy-2-(hydroxymethyl) tetrahydro-2H-pyran-3-yl)oxy)-3,5-dihydroxy-6-(hydroxymethyl) tetrahydro-2H-pyran-4-yl)oxy)-4,5-dihydroxy-6-(hydroxymethyl) tetrahydro-2H-pyran-3-yl) acetamide

Synonym: lyso-Globoside

MF: C₄₄H₈₀N₂O₂₂

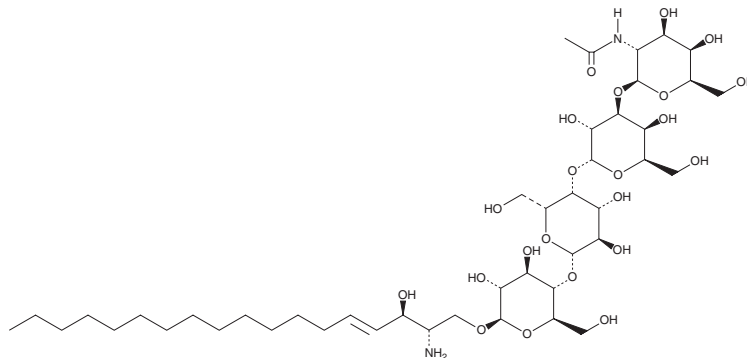
FW: 989.1

Purity: ≥98%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

Lyso-globotetraosylceramide (d18:1) is supplied as a solid. A stock solution may be made by dissolving the lyso-globotetraosylceramide (d18:1) in the solvent of choice, which should be purged with an inert gas. Lyso-globotetraosylceramide (d18:1) is soluble in a 2:1:0.1 solution of chloroform:methanol:water.

Description

Lyso-globotetraosylceramide is a form of globotetraosylceramide (Gb4) lacking the fatty acyl group. It binds to the Shiga-like toxin pig verotoxin of edema disease (VTE) in a glycolipid binding assay.¹ Lyso-globotetraosylceramide is acylated to stearic acid (Item No. 10011298) via a condensation reaction mediated by sphingolipid ceramide N-deacylase to form a ceramide.² As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. DeGrandis, S., Law, H., Brunton, J., et al. Globotetraosylceramide is recognized by the pig edema disease toxin. *J. Biol. Chem.* **264**(21), 12520-12525 (1989).
2. Kita, K., Kurita, T., and Ito, M. Characterization of the reversible nature of the reaction catalyzed by sphingolipid ceramide N-deacylase. A novel form of reverse hydrolysis reaction. *Eur. J. Biochem.* **268**(3), 592-602 (2001).

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, 10/30/2019

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