

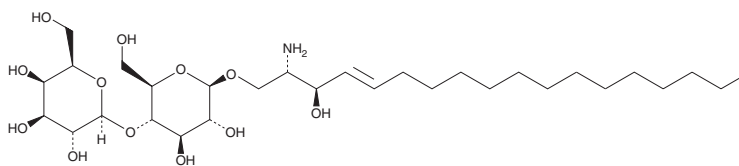
PRODUCT INFORMATION



Lactosylsphingosine (d18:1) (synthetic)

Item No. 24867

CAS Registry No.: 109785-20-8
Formal Name: (2S,3R,4E)-2-amino-3-hydroxy-4-octadecen-1-yl 4-O-β-D-galactopyranosyl-β-D-glucopyranoside
Synonym: Lactosyl Sphingosine, Lyso-Lactosylceramide (synthetic)
MF: C₃₀H₅₇NO₁₂
FW: 623.8
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

Lactosylsphingosine (d18:1) (synthetic) is supplied as a solid. A stock solution may be made by dissolving the Lactosylsphingosine (d18:1) (synthetic) in the solvent of choice, which should be purged with an inert gas. Lactosylsphingosine (d18:1) (synthetic) is soluble in DI water, DMSO, and a 2:1:0.1 solution of chloroform:methanol:DI water. We do not recommend storing the aqueous solution for more than one day.

Description

Lactosylsphingosine is a bioactive sphingolipid and a form of lactosylceramide (Item No. 16983) that is lacking the fatty acyl group. Lactosylsphingosine (1-50 μM) reduces viability of human neutrophils in a concentration-dependent manner.¹ Unlike lactosylceramide, lactosylsphingosine has no effect on protein synthesis and cell proliferation in cardiomyocytes.² Lactosylsphingosine is a precursor in the synthesis of lyso-ganglioside G_{M3}.³

References

1. Fiore, S., Nicolaou, K.C., Caulfield, T., *et al.* Evaluation of synthetic sphingosine, lysosphingolipids and glycosphingolipids as inhibitors of functional responses of human neutrophils. *Biochem. J.* **266**(1), 25-31 (1990).
2. Mishra, S. and Chatterjee, S. Lactosylceramide promotes hypertrophy through ROS generation and activation of ERK1/2 in cardiomyocytes. *Glycobiology* **24**(6), 518-531 (2014).
3. Thon, V., Lau, K., Yu, H., *et al.* PmST2: A novel *Pasteurella multocida* glycolipid α2-3-sialyltransferase. *Glycobiology* **21**(9), 1206-1216 (2011).

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, 11/15/2022

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