

PRODUCT INFORMATION



C22 Galactosylceramide (d18:1/22:0)

Item No. 9003458

CAS Registry No.: 74645-25-3
Formal Name: N-[(1S,2R,3E)-1-[(β-D-galactopyranosyloxy)methyl]-2-hydroxy-3-heptadecen-1-yl]-docosanamide

Synonyms: N-Docosanoyl-β-D-Galactosylceramide, Galactosylceramide (d18:1/22:0), GalCer(d18:1/22:0)

MF: C₄₆H₈₉NO₈

FW: 784.2

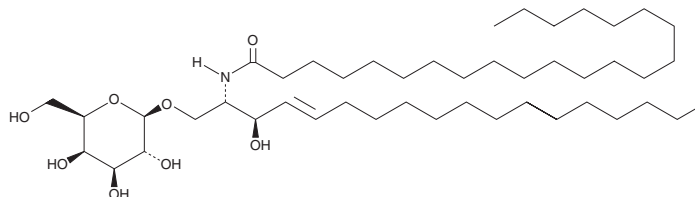
Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

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



Description

C22 Galactosylceramide is a sphingolipid that has been found in the central nervous system.^{1,2} C22 Galactosylceramide levels are increased in the spinal cord, but not the brain or liver, of mice fed a methionine-restricted diet.¹ C22 Galactosylceramide has been found in postmortem hippocampus from patients with Alzheimer's disease.² As this product is derived from a natural source, there may be variations in the sphingoid backbone.

References

1. Jové, M., Ayala, V., Ramírez-Núñez, O., *et al.* Specific lipidome signatures in central nervous system from methionine-restricted mice. *J. Proteome Res.* **12(6)**, 2679-2689 (2013).
2. Hejazi, L., Wong, J.W.H., Cheng, D., *et al.* Mass and relative elution time profiling: Two-dimensional analysis of sphingolipids in Alzheimer's disease brains. *Biochem. J.* **438(1)**, 165-175 (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, 10/28/2022

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