

PRODUCT INFORMATION



α C16 Galactosylceramide (d18:1/16:0)

Item No. 26009

CAS Registry No.: 2260795-77-3
Formal Name: N-[(1S,2R,3E)-1-[(α -D-galactopyranosyloxy)methyl]-2-hydroxy-3-heptadecen-1-yl]-hexadecanamide

Synonyms: Galactosylceramide (d18:1/16:0), α -GalCer(d18:1/16:0), D-galactosyl- α -1,1'-N-palmitoyl-D-erythro-Sphingosine

MF: C₄₀H₇₇NO₈

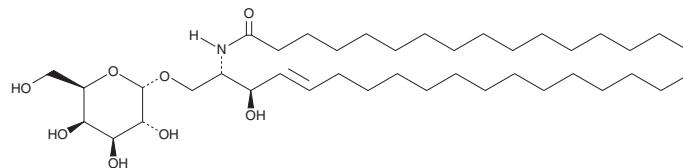
FW: 700.0

Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 2 years



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

Laboratory Procedures

α C16 Galactosylceramide (d18:1/16:0) is supplied as a solid. A stock solution may be made by dissolving the α C16 galactosylceramide (d18:1/16:0) in the solvent of choice, which should be purged with an inert gas. α C16 Galactosylceramide (d18:1/16:0) is slightly soluble in DMSO and methanol.

Description

α C16 Galactosylceramide is a glycosphingolipid that contains a galactose moiety attached to a ceramide acylated with palmitic acid (Item No. 10006627).

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, 10/06/2020

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