

PRODUCT INFORMATION



C18 Globotriaosylceramide-d₃ (d18:1/18:0-d₃)

Item No. 24626

Formal Name: N-((2S,3R,E)-1-(((2R,3R,4R,5S,6R)-5-(((2S,3R,4R,5R,6R)-3,4-dihydroxy-6-(hydroxymethyl)-5-(((2R,3R,4S,5R,6R)-3,4,5-trihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)tetrahydro-2H-pyran-2-yl)oxy)-3,4-dihydroxy-6-(hydroxymethyl)tetrahydro-2H-pyran-2-yl)oxy)-3-hydroxyoctadec-4-en-2-yl)octadecanamide-d₃

Synonyms: N-ω-CD₃-Octadecanoyl-ceramide Trihexoside, Galα1-4Galβ1-4Glcβ-Cer(d18:1/18:0) d₃, Gb3(d18:1/18:0-d₃), Globotriaosylceramide-d₃ (d18:1/18:0-d₃)

MF: C₅₄H₉₈D₃NO₁₈
FW: 1,055.4

Chemical Purity: ≥98% (C18 Globotriaosylceramide)

Deuterium

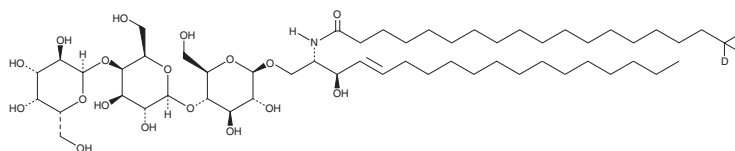
Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀

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

C18 Globotriaosylceramide-d₃ (d18:1/18:0-d₃) is supplied as a solid. A stock solution may be made by dissolving the C18 globotriaosylceramide-d₃ (d18:1/18:0-d₃) in the solvent of choice. C18 Globotriaosylceramide-d₃ (d18:1/18:0-d₃) is soluble in DMSO and in a 2:1 solution of chloroform:methanol.

Description

C18 Globotriaosylceramide-d₃ is intended for use as an internal standard for the quantification of C18 globotriaosylceramide by GC- or LC-MS. C18 globotriaosylceramide is an endogenous sphingolipid that accumulates in the brain, heart, kidney, liver, lung, and spleen in a mouse model of Fabry disease, a lysosomal storage disorder characterized by a deficiency in the enzyme α-galactosidase A.¹ As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Reference

1. Kamani, M.A., Provençal, P., Boutin, M., et al. Glycosphingolipid storage in Fabry mice extends beyond globotriaosylceramide and is affected by ABCB1 depletion. *Future Sci. OA.* **2(4):FS0147**, (2016).

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, 11/18/2022

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