

PRODUCT INFORMATION



C23 Globotriaosylceramide (d18:1/23:0)

Item No. 24878

CAS Registry No.: 536745-84-3

Formal Name: N-[(1S,2R,3E)-1-[[[O- α -D-galactopyranosyl-(1 \rightarrow 4)-O- β -D-galactopyranosyl-(1 \rightarrow 4)- β -D-glucopyranosyl]oxy]methyl]-2-hydroxy-3-heptadecen-1-yl]-tricosanamide

Synonyms: C23 Ceramide Trihexoside (d18:0/23:0), N-Tricosanoyl-ceramide Trihexoside, N-Tricosanoyl Globotriaosylceramide

MF: C₅₉H₁₁₁NO₁₈

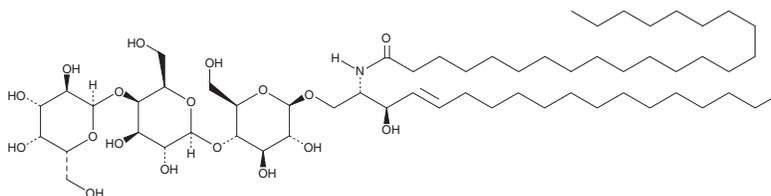
FW: 1,122.5

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



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

Laboratory Procedures

C23 Globotriaosylceramide (d18:1/23:0) is supplied as a solid. A stock solution may be made by dissolving the C23 globotriaosylceramide (d18:1/23:0) in the solvent of choice, which should be purged with an inert gas. C23 Globotriaosylceramide (d18:1/23:0) is soluble in DMSO, methanol (hot), and a 2:1 solution of chloroform:methanol.

Description

C23 Globotriaosylceramide is a sphingolipid that has been used as a standard for the quantification of globotriaosylceramide (Item Nos. 24870 | 24871 | 24872) in human fibroblasts by HPLC/APCI-MS.¹ As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Reference

1. Farwanah, H., Wirtz, J., Kolter, T., et al. Normal phase liquid chromatography coupled to quadrupole time of flight atmospheric pressure chemical ionization mass spectrometry for separation, detection and mass spectrometric profiling of neutral sphingolipids and cholesterol. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* **877**(27), 2976-2982 (2009).

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, 10/27/2022

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