

PRODUCT INFORMATION



N-Glycine Globotriaosylsphingosine (d18:1)

Item No. 24874

CAS Registry No.: 1360882-59-2

Formal Name: 2-amino-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]-acetamide

Synonyms: N-Glycinated Globotriaosylsphingosine, N-Glycinated Lyso-Ceramide Trihexoside, N-Glycine Lyso-Globotriaosylceramide (d18:1)

MF: $C_{38}H_{70}N_2O_{18}$

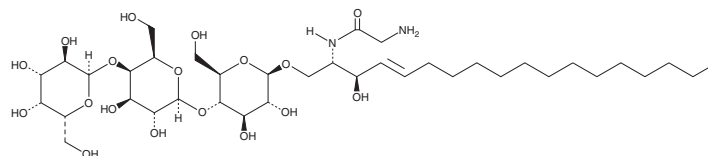
FW: 843.0

Purity: $\geq 98\%$

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

N-Glycine Globotriaosylsphingosine (d18:1) is supplied as a solid. A stock solution may be made by dissolving the N-Glycine Globotriaosylsphingosine (d18:1) in the solvent of choice. N-Glycine Globotriaosylsphingosine (d18:1) is soluble in dimethyl formamide and a 2:1:0.1 solution of chloroform:methanol:DI water.

Description

N-Glycine Globotriaosylsphingosine is a glycinated form of globotriaosylceramide (Item Nos. 24870 | 24871 | 24872) that contains a glycine residue in place of the fatty acyl group. It has been used as an internal standard for the quantification of lyso-globotriaosylceramide (Item No. 24873) in human plasma by UPLC-MS/MS.¹ As this product is derived from a natural source, there may be variations in the sphingoid backbone.

Reference

1. Krüger, R., Tholey, A., Jakoby, T., et al. Quantification of the Fabry marker lysoGb3 in human plasma by tandem mass spectrometry. *J. Chromatogr. B. Analyt. Technol. Biomed. Life Sci.* **883-884**, 128-135 (2012).

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, 10/31/2022

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