

# PRODUCT INFORMATION



## C17 Globotriaosylceramide (d18:1/17:0)

Item No. 24876

CAS Registry No.: 536745-81-0

Formal Name: N-[(1S,2R,3E)-1-[[[O-α-D-galactopyranosyl-(1→4)-O-β-D-galactopyranosyl-(1→4)-β-D-glucopyranosyl]oxy)methyl]-2-hydroxy-3-heptadecenyl]-heptadecanamide

Synonyms: C17 Ceramide Trihexoside (d18:1/17:0), C17:0 Gb<sub>3</sub>, N-Heptadecanoyl Ceramide Trihexoside

MF: C<sub>53</sub>H<sub>99</sub>NO<sub>18</sub>

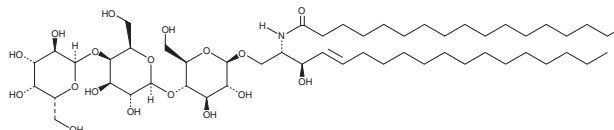
FW: 1,038.3

Purity: ≥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

C17 Globotriaosylceramide (d18:1/17:0) is supplied as a solid. A stock solution may be made by dissolving the C17 globotriaosylceramide (d18:1/17:0) in the solvent of choice, which should be purged with an inert gas. C17 Globotriaosylceramide (d18:1/17:0) is soluble in DMSO, methanol (hot), and a 2:1 solution of chloroform:methanol.

### Description

C17 Globotriaosylceramide is a sphingolipid that has been used as an internal standard for the quantification of globotriaosylceramide (Item Nos. 24870 | 24871 | 24872) in plasma and urine from patients with Fabry disease by GC-MS.<sup>1</sup> As this product is derived from a natural source, there may be variations in the sphingoid backbone.

### Reference

1. Winchester, B. and Mills, K. Internal standards for sphingolipids for use in mass spectrometry. *Ich Productions Limited WO03/048784A2* (2001).

#### 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, 07/14/2023

#### 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