

# PRODUCT INFORMATION



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

Item No. 24356

**CAS Registry No.:** 1354699-26-5  
**Formal Name:** N-[(1S,2R,3E)-1-[[[4-O-β-D-galactopyranosyl]-β-D-glucopyranosyl]oxy]methyl]-2-hydroxy-3-heptadecen-1-yl]-heptadecanamide

**Synonyms:** N-Heptadecanoyl Lactosylceramide, LacCer(d18:1/17:0), Lactosylceramide (d18:1/17:0)

**MF:** C<sub>47</sub>H<sub>89</sub>NO<sub>13</sub>

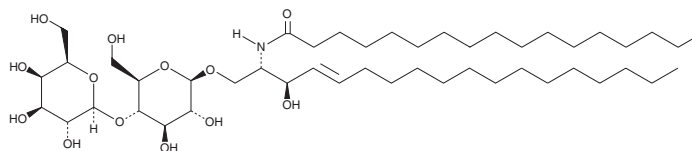
**FW:** 876.2

**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 Lactosylceramide (d18:1/17:0) is supplied as a solid. A stock solution may be made by dissolving the C17 lactosylceramide (d18:1/17:0) in the solvent of choice. C17 Lactosylceramide (d18:1/17:0) is soluble in a 2:1:0.1 solution of chloroform:methanol:water.

### Description

C17 Lactosylceramide is a naturally occurring sphingolipid that has been found in human brain metastases of lung adenocarcinomas but not in healthy brain tissue.<sup>1</sup> As this product is derived from a natural source, there may be variations in the sphingoid backbone.

### Reference

1. Zamfir, A.D., Serb, A., Vukeli, Ž., *et al.* Assessment of the molecular expression and structure of gangliosides in brain metastasis of lung adenocarcinoma by an advanced approach based on fully automated chip-nanoelectrospray mass spectrometry. *J. Am. Soc. Mass Spectrom.* **22(12)**, 2145-2159 (2011).

#### 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, 12/06/2022

#### 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