

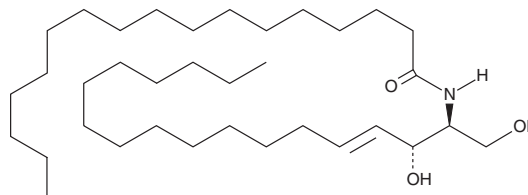
PRODUCT INFORMATION



C17 Ceramide (d18:1/17:0)

Item No. 22532

CAS Registry No.: 67492-16-4
Formal Name: N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-heptadecanamide
Synonyms: Cer(d18:1/17:0), Ceramide (d18:1/17:0), N-heptadecanoyl-D-erythro-Sphingosine
MF: C₃₅H₆₉NO₃
FW: 551.9
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 Ceramide (d18:1/17:0) is supplied as a solid. A stock solution may be made by dissolving the C17 ceramide (d18:1/17:0) in the solvent of choice, which should be purged with an inert gas. C17 Ceramide (d18:1/17:0) is soluble in organic solvents such as ethanol (heated), methanol (heated), and chloroform.

Description

C17 Ceramide is a synthetic ceramide that contains a non-natural C17:0 fatty acid acylated to sphingosine.¹ It has been used as a marker of arylsulfatase A and sphingolipid activator B activity in patients undergoing enzyme replacement therapy for the genetic disease metachromatic leukodystrophy.

Reference

1. Cui, Y., Colsch, B., Afonso, C., *et al.* Synthetic sulfogalactosylceramide (sulfatide) and its use for the mass spectrometric quantitative urinary determination in metachromatic leukodystrophies. *Glycoconj. J.* **25(2)**, 147-155 (2008).

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, 12/02/2022

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