

PRODUCT INFORMATION



C16 Ceramide (d18:1(14Z)/16:0)

Item No. 28601

Formal Name: N-((2S,3R,Z)-1,3-dihydroxyoctadec-14-en-2-yl)palmitamide

Synonyms: C16:0-Cer, Cer(d18:1(14Z)/16:0), Ceramide (d18:1(14Z)/16:0), N-palmitoyl-4,5-dihydro-14Z-Sphingosine

MF: C₃₄H₆₇NO₃

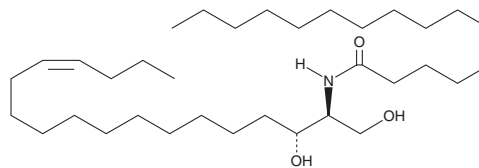
FW: 537.9

Purity: ≥95%

Supplied as: A crystalline 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

C16 Ceramide (d18:1(14Z)/16:0) is supplied as a crystalline solid. A stock solution may be made by dissolving the C16 ceramide (d18:1(14Z)/16:0) in the solvent of choice, which should be purged with an inert gas. C16 Ceramide (d18:1(14Z)/16:0) is soluble in the organic solvent dimethyl formamide at a concentration of approximately 0.5 mg/ml.

Description

C16 Ceramide (d18:1(14Z)/16:0) is a long-chain atypical ceramide containing a sphingosine (d18:1(14Z)) (Item No. 27019) backbone with a *cis* double bond at the 14-15 position rather than a *trans* double bond at the 4-5 position.

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, 03/19/2024

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