

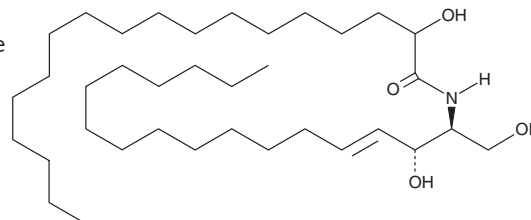
PRODUCT INFORMATION



C18 ((±)-2'-hydroxy) Ceramide (d18:1/18:0)

Item No. 22687

CAS Registry No.: 34249-41-7
Formal Name: 2-hydroxy-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-octadecanamide
Synonym: ((±)-2'-hydroxy) Cer(d18:1/18:0), ((±)-2'-hydroxy) Ceramide (d18:1/18:0), 18:0 (2R/S-OH)-Ceramide,N-(2'-(R,S)-hydroxystearoyl)-D-erythro-Sphingosine
MF: C₃₆H₇₁NO₄
FW: 582.0
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

C18 ((±)-2'-hydroxy) Ceramide (d18:1/18:0 (18:0 (2R/S-OH)-ceramide) is supplied as a solid. A stock solution may be made by dissolving the 18:0 (2R/S-OH)-ceramide in the solvent of choice. 18:0 (2R/S-OH)-Ceramide is soluble in a 4:3:1 solution of chloroform:methanol:water.

Description

18:0 (2R/S-OH)-Ceramide is a bioactive sphingolipid found in the stratum corneum layer of mammalian epidermis.¹ 18:0 (2R/S-OH)-Ceramide plays an important role in maintaining normal hydration levels for physiological function of the epidermis.

Reference

1. Tfayli, A., Jamal, D., Vyumvuhore, R., *et al.* Hydration effects on the barrier function of stratum corneum lipids: Raman analysis of ceramides 2, III and 5. *Analyst* **138(21)**, 6582-6588 (2013).

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, 11/14/2022

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