

PRODUCT INFORMATION



C12((±)-2'-hydroxy) dihydro Ceramide (d18:0/12:0)

Item No. 24435

CAS Registry No.: 2377379-54-7

Formal Name: 2-hydroxy-N-[(1S,2R)-2-hydroxy-1-(hydroxymethyl)heptadecyl]-dodecanamide

Synonyms: (±)-2'-hydroxy Cer(d18:0/12:0), (±)-2'-hydroxy Ceramide (d18:0/12:0), N-(R,S)-α-hydroxydodecanoyl-D-erythro-Dihydrosphingosine

MF: C₃₀H₆₁NO₄

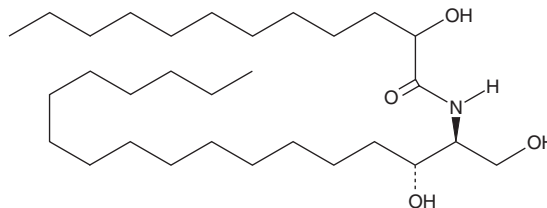
FW: 499.8

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

C12((±)-2'-hydroxy) dihydro Ceramide (d18:0/12:0) is supplied as a solid. A stock solution may be made by dissolving the C12((±)-2'-hydroxy) dihydro ceramide (d18:0/12:0) in the solvent of choice, which should be purged with an inert gas. C12((±)-2'-hydroxy) dihydro Ceramide (d18:0/12:0) is soluble in a 2:1:0.5 solution of chloroform:methanol:water.

Description

C12((±)-2'-hydroxy) dihydro Ceramide (d18:0/12:0) is a 2'-hydroxylated form of C12 dihydro ceramide.

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, 01/06/2023

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