

PRODUCT INFORMATION



C24 (2'(R)-hydroxy) Ceramide (d18:1/24:0)

Item No. 22527

CAS Registry No.: 64655-47-6
Formal Name: (2R)-2-hydroxy-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-tetracosanamide
Synonyms: 2'(R)-hydroxy Cer(d18:1/24:0),
2'-hydroxy C24-Ceramide,
(2'(R)-hydroxy) Ceramide (d18:1/24:0),
N-(2'(R)-hydroxylignoceroyl)-D-erythro-Sphingosine,
N-(R)- α -Hydroxytetracosanoyl-D-erythro-sphingosine

MF: C₄₂H₈₃NO₄

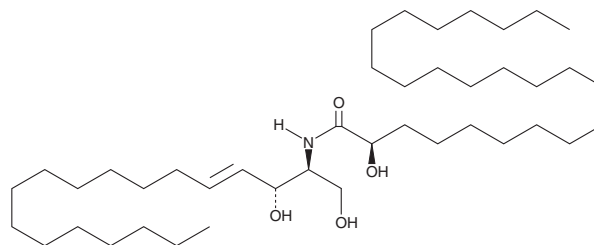
FW: 666.1

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

C24 (2'(R)-hydroxy) Ceramide (d18:1/24:0) is supplied as a solid. A stock solution may be made by dissolving the C24 (2'(R)-hydroxy) ceramide (d18:1/24:0) in the solvent of choice, which should be purged with an inert gas. C24 (2'(R)-hydroxy) Ceramide (d18:1/24:0) is soluble in a 4:1 solution of chloroform:methanol. C24 (2'(R)-hydroxy) Ceramide (d18:1/24:0) is also soluble in ethanol (warmed).

Description

C24 (2'(R)-hydroxy) Ceramide is a synthetic ceramide and an isomer of C24 (2'(S)-hydroxy) ceramide (Item No. 22526).¹ It has been used in the preparation of synthetic lipid membranes that model lamellar and lateral stratum corneum lipid organization.

Reference

1. de Jager, M.W., Gooris, G.S., Ponec, M., *et al.* Lipid mixtures prepared with well-defined synthetic ceramides closely mimic the unique stratum corneum lipid phase behavior. *J. Lipid Res.* **46**(12), 2649-2656 (2005).

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, 11/18/2022

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