

# PRODUCT INFORMATION



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

Item No. 28935

**CAS Registry No.:** 164989-36-0  
**Formal Name:** (2R)-2-hydroxy-N-[(1S,2R)-2-hydroxy-1-(hydroxymethyl)heptadecyl]-tetracosanamide  
**Synonyms:** (R)-2'-hydroxy Cer(d18:0/24:0), (R)-2'-hydroxy Ceramide (d18:0/24:0), N-(R)- $\alpha$ -C24 dihydro Ceramide (d18:0/24:0), N-(R)- $\alpha$ -hydroxytetracosanoyl-D-erythro-Dihydrosphingosine

**MF:** C<sub>42</sub>H<sub>85</sub>NO<sub>4</sub>

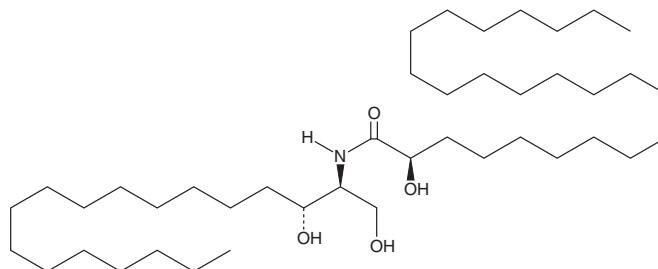
**FW:** 668.1

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

### Description

C24 (2'(R)-hydroxy) dihydro ceramide is the (2R) isomer of a 2'-hydroxylated form of C24 dihydro ceramide (Item No. 25602) and the (R) isomer of C24 (2'(S)-hydroxy) dihydro ceramide (Item No. 28936).

#### 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, 12/06/2022

#### 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