

# PRODUCT INFORMATION



## C16 (2'(S)-hydroxy) Ceramide (d18:1/16:0)

Item No. 22516

**CAS Registry No.:** 890041-50-6  
**Formal Name:** (2S)-2-hydroxy-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-hexadecanamide

**Synonyms:** 16:0 (2S-OH)-Ceramide, 2'-hydroxy C16-Ceramide

**MF:** C<sub>34</sub>H<sub>67</sub>NO<sub>4</sub>

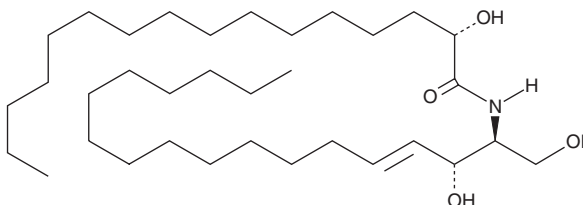
**FW:** 553.9

**Purity:** ≥95%

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥2 years



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

### Laboratory Procedures

C16 (2'(S)-hydroxy) Ceramide (d18:1/16:0) is supplied as a crystalline solid. A stock solution may be made by dissolving the C16 (2'(S)-hydroxy) ceramide (d18:1/16:0) in the solvent of choice. C16 (2'(S)-hydroxy) Ceramide (d18:1/16:0) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. C16 (2'(S)-hydroxy) Ceramide (d18:1/16:0) is miscible in ethanol and has a solubility of approximately 2 and 10 mg/ml in DMSO and DMF, respectively.

### Description

C16 (2'(S)-hydroxy) Ceramide is the synthetic ceramide stereoisomer of N-(2'(R)-hydroxypalmitoyl)-D-erythro-sphingosine.<sup>1</sup> C16 (2'(S)-hydroxy) Ceramide is the less active isomer, having no effect on C6 glioma cell viability at concentrations up to 20 μM.

### Reference

1. Kota, V., Dhople, V.M., Fullbright, G., *et al.* 2'-hydroxy C16-ceramide induces apoptosis-associated proteomic changes in C6 glioma cells. *J. Proteome Res.* **12(10)**, 4366-4375 (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, 01/25/2018

#### 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