

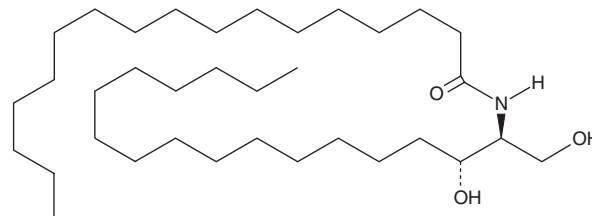
# PRODUCT INFORMATION



## C17 dihydro Ceramide (d18:0/17:0)

Item No. 24427

**CAS Registry No.:** 1388156-40-8  
**Formal Name:** N-[(1S,2R)-2-hydroxy-1-(hydroxymethyl)heptadecyl]-heptadecanamide  
**Synonyms:** Cer(d18:0/17:0), Ceramide (d18:0/17:0), N-heptadecanoyl-D-erythro-Dihydrosphingosine  
**MF:** C<sub>35</sub>H<sub>71</sub>NO<sub>3</sub>  
**FW:** 553.9  
**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

C17 dihydro Ceramide (d18:0/17:0) is supplied as a solid. A stock solution may be made by dissolving the C17 dihydro ceramide (d18:0/17:0) in the solvent of choice. C17 dihydro Ceramide (d18:0/17:0) is soluble in ethanol (heated) and DMSO.

### Description

C17 dihydro Ceramide is a synthetic ceramide that contains a non-natural C17:0 fatty acid acylated to sphingosine.<sup>1,2</sup> Conversion of C17 dihydro ceramide to C17 ceramide (Item No. 22532) has been used to study dihydroceramide desaturase activity and ceramide metabolism in cells.

### References

1. Spassieva, S., Seo, J.G., Jiang, J.C., *et al.* Necessary role for the Lag1p motif in (dihydro)ceramide synthase activity. *J. Biol. Chem.* **281**(45), 33931-33938 (2006).
2. Spassieva, S.D., Rahmaniyan, M., Bielawski, J., *et al.* Cell density-dependent reduction of dihydroceramide desaturase activity in neuroblastoma cells. *J. Lipid Res.* **53**(5), 918-928 (2012).

#### 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, 11/29/2022

#### 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