

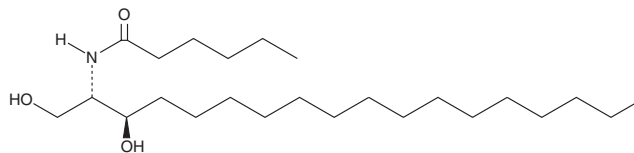
# PRODUCT INFORMATION



## C6 dihydro Ceramide (d18:0/6:0)

Item No. 24368

**CAS Registry No.:** 171039-13-7  
**Formal Name:** N-[(1S,2R)-2-hydroxy-1-(hydroxymethyl)heptadecyl]-hexanamide  
**Synonyms:** Cer(d18:0/6:0), Ceramide (d18:0/6:0), N-hexanoyl-D-erythro-Dihydrosphingosine  
**MF:** C<sub>24</sub>H<sub>49</sub>NO<sub>3</sub>  
**FW:** 399.7  
**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

C6 dihydro Ceramide (d18:0/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 dihydro ceramide (d18:0/6:0) in the solvent of choice. C6 dihydro Ceramide (d18:0/6:0) is soluble in organic solvents such as chloroform, ethanol, and DMSO, which should be purged with an inert gas.

### Description

C6 dihydro Ceramide is a metabolically inactive analog of C6 ceramide (Item No. 62525) that has been used as a negative control in experiments investigating the biological activity of C6 ceramide.<sup>1</sup>

### Reference

1. Cai, Z., Bettaieb, A., El Mahdani, N., *et al.* Alteration of the sphingomyelin/ceramide pathway is associated with resistance of human breast carcinoma MCF7 cells to tumor necrosis factor- $\alpha$ -mediated cytotoxicity. *J. Biol. Chem.* **272**(11), 6918-6926 (1997).

**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, 12/19/2022

### 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](http://WWW.CAYMANCHEM.COM)