

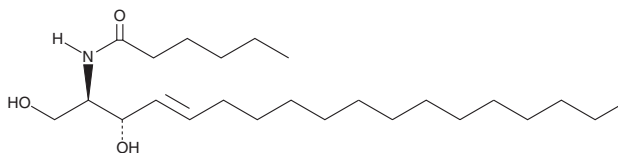
# PRODUCT INFORMATION



## C6 L-erythro Ceramide (d18:1/6:0)

Item No. 24388

**CAS Registry No.:** 189894-78-8  
**Formal Name:** N-[(1R,2S,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-hexanamide  
**Synonyms:** L-erythro Cer(d18:1/6:0), L-erythro Ceramide (d18:1/6:0), N-hexanoyl-L-erythro-Sphingosine  
**MF:** C<sub>24</sub>H<sub>47</sub>NO<sub>3</sub>  
**FW:** 397.6  
**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 L-erythro Ceramide (d18:1/6:0) is supplied as a solid. A stock solution may be made by dissolving the C6 L-erythro ceramide (d18:1/6:0) in the solvent of choice. C6 L-erythro Ceramide (d18:1/6:0) is soluble in organic solvents such as chloroform, ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of C6 L-erythro ceramide (d18:1/6:0) in DMF is approximately 5 mg/ml.

### Description

C6 L-erythro Ceramide is a bioactive sphingolipid and cell-permeable analog of naturally occurring ceramides.<sup>1,2</sup> It is metabolized by ceramide glucosyltransferase to form C6 L-erythro glucosylceramide. C6 L-erythro Ceramide is cytotoxic to U937 cells (IC<sub>50</sub> = 18 μM).<sup>3</sup>

### References

- Schwarz, A. and Futerman, A.H. Distinct roles for ceramide and glucosylceramide at different stages of neuronal growth. *J. Neurosci.* **17**(9), 2929-2938 (1997).
- Paul, P., Kamisaka, Y., Marks, D.L., et al. Purification and characterization of UDP-glucose: Ceramide glucosyltransferase from rat liver Golgi membranes. *J. Biol. Chem.* **271**(4), 2287-2293 (1996).
- Chang, Y.-T., Choi, J., Ding, S., et al. The synthesis and biological characterization of a ceramide library. *J. Am. Chem. Soc.* **124**(9), 1856-1857 (2002).

#### 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, 10/31/2022

#### 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