

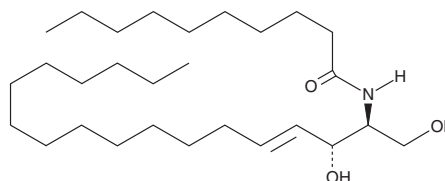
# PRODUCT INFORMATION



## C10 Ceramide (d18:1/10:0)

Item No. 22529

**CAS Registry No.:** 111122-57-7  
**Formal Name:** N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-decanamide  
**Synonyms:** Cer(d18:1/10:0), N-decanoyl-D-erythro-Sphingosine, C10 Ceramide, Ceramide (d18:1/10:0), N-decanoyl-D-erythro-Sphingosine  
**MF:** C<sub>28</sub>H<sub>55</sub>NO<sub>3</sub>  
**FW:** 453.7  
**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

C10 Ceramide (d18:1/10:0) is supplied as a solid. A stock solution may be made by dissolving the C10 ceramide (d18:1/10:0) in the solvent of choice, which should be purged with an inert gas. C10 Ceramide (d18:1/10:0) is soluble in the organic solvent dimethyl formamide at a concentration of approximately 0.50 mg/ml.

### Description

C10 Ceramide is an endogenous, bioactive sphingolipid produced by ceramide synthase.<sup>1</sup> C10 ceramide inhibits growth of wild-type and MDR1 gene-transfected MDA435/LCC6 human breast cancer cells with IC<sub>50</sub> values of 44.1 and 46 μM, respectively. It also inhibits the growth of J774 murine macrophages in a dose-dependent manner.

### Reference

1. Shabbits, J.A. and Mayer, L.D. Intracellular delivery of ceramide lipids via liposomes enhances apoptosis in vitro. *Biochim Biophys. Acta.* **1612(1)**, 98-106 (2003).

#### 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, 05/10/2023

#### 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