

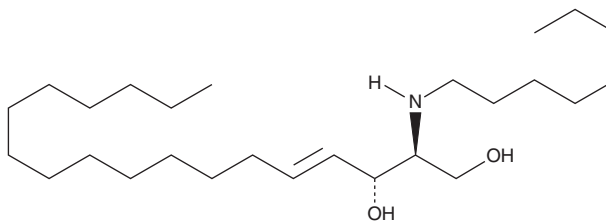
# PRODUCT INFORMATION



## C8 Ceramine (d18:1/8:0)

Item No. 62545

**CAS Registry No.:** 170926-06-4  
**Formal Name:** N-octylsphingosine  
**Synonyms:** C8 Ceramine, Ceramine (d18:1/8:0)  
**MF:** C<sub>26</sub>H<sub>53</sub>NO<sub>2</sub>  
**FW:** 411.7  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

C8 Ceramine is supplied as a crystalline solid. A stock solution may be made by dissolving the C8 ceramine in the solvent of choice. C8 Ceramine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of C8 ceramine in these solvents is approximately 33, 20, and 22 mg/ml, respectively.

C8 Ceramine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, C8 ceramine should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. C8 Ceramine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

C8 Ceramine is an analog of C8 ceramide (Item No. 62540) in which the carbonyl group of the ceramide is replaced by a methylene group. At a concentration of 10 μM, C8 ceramine induces maximal DNA fragmentation in U937 cells after 6 hours of incubation, compared to 12 hours for C8 ceramide.<sup>1</sup>

### Reference

1. Bernert, J.T., Jr., Bell, C.J., McGuffey, J.E., *et al.* Determination of "free" glycerol in human serum reference materials by isotope-dilution gas chromatography-mass spectrometry. *Journal of Chromatography* **578(1)**, 1-7 (1992).

**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, 02/20/2018

### 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