

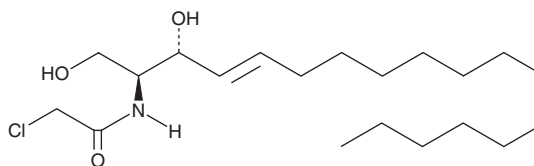
# PRODUCT INFORMATION



## SOCLAC

Item No. 9003928

**CAS Registry No.:** 126586-82-1  
**Formal Name:** 2-chloro-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-acetamide  
**Synonyms:** N-chloroacetyl Sphingosine, PKS2  
**MF:** C<sub>20</sub>H<sub>38</sub>ClNO<sub>3</sub>  
**FW:** 376.0  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

SOCLAC is supplied as a solid. A stock solution may be made by dissolving the SOCLAC in the solvent of choice, which should be purged with an inert gas. SOCLAC is soluble in the organic solvent chloroform.

### Description

SOCLAC is an irreversible inhibitor of acid ceramidase ( $K_i = 40.2 \text{ nM}$ ) and a derivative of the acid ceramidase inhibitor SABRAC (Item No. 9003925).<sup>1</sup>

### Reference

1. Ordóñez, Y.F., Abad, J.L., Aseeri, M., *et al.* Activity-based imaging of acid ceramidase in living cells. *J. Am. Chem. Soc.* **141**(19), 7736-7742 (2019).

#### 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, 01/06/2022

#### 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)