

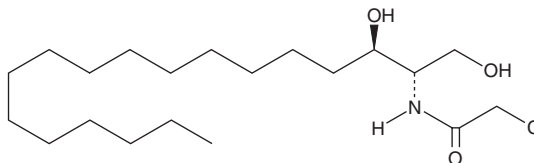
PRODUCT INFORMATION



SACLAC

Item No. 9003926

CAS Registry No.: 2248703-42-4
Formal Name: 2-chloro-N-[(1S,2R)-2-hydroxy-1-(hydroxymethyl)heptadecyl]-acetamide
MF: $C_{20}H_{40}ClNO_3$
FW: 378.0
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

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

Description

SACLAC is an irreversible inhibitor of acid ceramidase ($K_i = 97.1$ nM) and a derivative of the acid ceramidase inhibitor SABRAC (Item No. 9003925).¹ It decreases levels of sphingosine-1-phosphate (S1P) and increases total ceramide levels in OCI-AML-2 acute myeloid leukemia (AML) cells when used at a concentration of $2.5 \mu M$.² SACLAC (10 and $20 \mu M$) induces apoptosis in primary AML cells. It reduces leukemic burden in MV4-11 and U937 AML mouse xenograft models when administered at a dose of 5 mg/kg.

References

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).
2. Pearson, J.M., Tan, S.-F., Sharma, A., *et al.* Ceramide analogue SACLAC modulates sphingolipid levels and MCL-1 splicing to induce apoptosis in acute myeloid leukemia. *Mol. Cancer Res.* **18**(3), 352-363 (2020).

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