

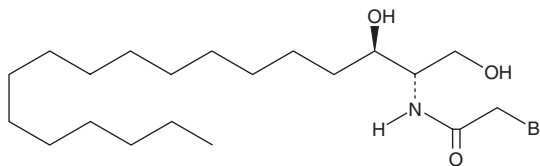
PRODUCT INFORMATION



SABRAC

Item No. 9003925

CAS Registry No.: 886212-98-2
Formal Name: 2-bromo-N-[(1S,2R)-2-hydroxy-1-(hydroxymethyl)heptadecyl]-acetamide
MF: $C_{20}H_{40}BrNO_3$
FW: 422.4
Purity: $\geq 95\%$
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

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

Description

SABRAC is an inhibitor of acid ceramidase ($IC_{50} = 52 \text{ nM}$).¹ It is selective for acid ceramidase over neutral ceramidase and the cysteine protease papain at $3 \mu\text{M}$. SABRAC inhibits acid ceramidase in PC3 MC metastatic prostate cancer cells ($IC_{50} = <1 \mu\text{M}$), as well as induces accumulation of ceramides in the same cells. It completely inhibits anchorage-independent colony formation but does not induce cytotoxicity in PC3 MC cells when used at a concentration of $5 \mu\text{M}$.

Reference

1. Camacho, L., Meca-Cortés, O., Abad, J.L., *et al.* Acid ceramidase as a therapeutic target in metastatic prostate cancer. *J. Lipid Res.* **54**(5), 1207-1220 (2013).

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, 11/14/2025

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