

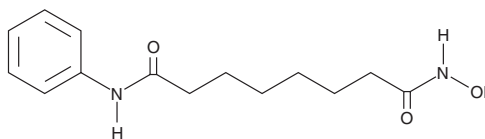
PRODUCT INFORMATION



SAHA

Item No. 10009929

CAS Registry No.: 149647-78-9
Formal Name: N¹-hydroxy-N⁸-phenyl-octanediamide
Synonyms: Suberoylanilide Hydroxamic Acid, Vorinostat
MF: C₁₄H₂₀N₂O₃
FW: 264.3
Purity: ≥98%
UV/Vis.: λ_{max}: 243 nm
Supplied as: A crystalline 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

SAHA is supplied as a crystalline solid. A stock solution may be made by dissolving the SAHA in an organic solvent purged with an inert gas. SAHA is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SAHA in ethanol is approximately 0.25 mg/ml and approximately 20 mg/ml in DMSO and DMF.

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

Description

SAHA is a histone deacetylase (HDAC) inhibitor that binds directly to the catalytic site of the enzyme thereby blocking substrate access. It inhibits class I, II, and IV HDACs at 50-200 nM and arrests cell growth in a wide variety of transformed cells in culture at 2.5-5.0 μM.¹

Reference

1. Marks, P.A. and Breslow, R. Dimethyl sulfoxide to vorinostat: Development of this histone deacetylase inhibitor as an anticancer drug. *Nature Biotechnology* **25(1)**, 84-90 (2007).

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, 08//17/2017

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