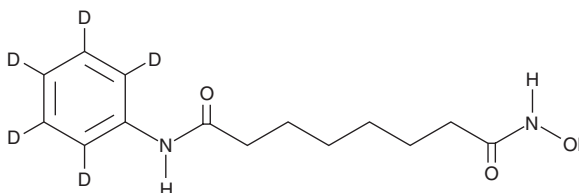


PRODUCT INFORMATION



SAHA-d₅
Item No. 22107

CAS Registry No.: 1132749-48-4
Formal Name: N¹-hydroxy-N⁸-(phenyl-2,3,4,5,6-d₅)-octanediamide
Synonyms: Suberoylanilide Hydroxamic Acid-d₅, Vorinostat-d₅
MF: C₁₄H₁₅D₅N₂O₃
FW: 269.4
Chemical Purity: ≥96% (SAHA)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₅); ≤1% d₀
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

SAHA-d₅ contains five deuterium atoms at the 2, 3, 4, 5, and 6 positions. It is intended for use as an internal standard for the quantification of SAHA (suberoylanilide hydroxamic acid; Item No. 10009929) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

SAHA-d₅ is supplied as a solid. A stock solution may be made by dissolving the SAHA-d₅ in the solvent of choice. SAHA-d₅ is slightly soluble in acetonitrile and DMSO.

Description

SAHA is a histone deacetylase (HDAC) inhibitor that binds directly to the catalytic site of the enzyme thereby blocking substrate access.¹ 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.

Reference

1. Marks, P.A. and Breslow, R. Dimethyl sulfoxide to vorinostat: Development of this histone deacetylase inhibitor as an anticancer drug. *Nat. Biotechnol.* **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, 12/06/2022

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