
SAHA

(Catalog No. M41000)

Alternative Names

Vorinostat, Suberanilohydroxamic Acid, MK-0683

Description

Potent, reversible inhibitor of histone deacetylase 1 (HDAC1) and 3 (HDAC3). Induces cell growth arrest at both G1 and G2 phases. Induces apoptosis.

CAS Number

149647-78-9

Chemical Formula

$C_{14}H_{20}N_2O_3$

Molecular Weight

264.3

Size

10 mg

Purity

≥99%

Solubility

DMSO (60 mg/ml) of EtOH (2 mg/ml)

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
SAHA	10 mg	M41000 -2

This product is for research purposes only. Not intended for use in diagnostic procedures.