

SAHA (Vorinostat)

Catalog Number P004-50MG

Catalog Number P004-250MG



ARBOR
ASSAYS

FEATURES

HDAC inhibitor

Cutaneous T-cell lymphoma treatment

In vitro and *in vivo* effects on latent HIV infected T-Cells



INTRODUCTION

SAHA (suberoylanilide hydroxamic acid, Vorinostat, Zolinza™) is a histone deacetylase (HDAC) inhibitor that binds directly to the catalytic site of the enzyme thereby blocking substrate access. Sirt activator based on a dihydropyridine structural scaffold. EC150 (effective concentration able to increase enzyme activity by 150%) is 1 μM for Sirt1, 25 μM for Sirt2, and 50 μM for Sirt3. The compound is active in whole cells and arrests cell cycle at G1/S phase in U937 cells. SAHA inhibits class I and class II HDACs at around 50 nM and arrests cell growth in a wide variety of transformed cells in culture at 2.5-5.0 μM.

FORM: White Powder

MOLECULAR WEIGHT: 264.3

STORAGE: Room temperature, desiccated for up to 1 year. Store solutions at -20°C for up to 6 months.

FORMULA: C₁₄H₂₀N₂O₃

CAS NUMBER: 149647-78-9

OTHER NAMES: N¹-hydroxy-N⁸-phenyl-octanediamide, suberoylanilide hydroxamic acid, Vorinostat, Zolinza™

USES: Soluble to 50 mg/mL in DMSO and 2 mg/mL in Ethanol.

REFERENCES:

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).