

## SAHA (Vorinostat)

**Catalog #:** 27006

**Size:** 100 mg

**CAS Registry #:** 149647-78-9

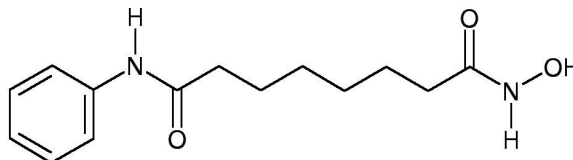
**Purity:** ≥ 95%

**Chemical Formula:** C<sub>14</sub>H<sub>20</sub>N<sub>2</sub>O<sub>3</sub>

**Molecular Weight:** 264.32

**Lot #:** 111118

**Structure:**



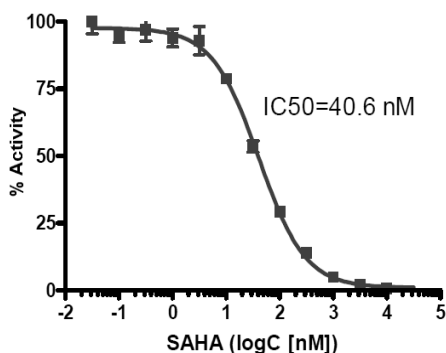
**Description:** Soruberoylanilide Hydroxamic Acid (SAHA), or Vorinostat, is a histone deacetylase inhibitor and causes growth arrest and death of some transformed cells both *in vitro* and *in vivo*, with little or no toxic effects on normal cells.

**Appearance:** White to off-white crystalline solid; granular or powder.

**Solubility:** Soluble in DMSO at 66 mg/mL; soluble in ethanol at 2 mg/mL with slight warming; very poorly soluble in water; maximum solubility in plain water is estimated to be about 20-50 μM; buffers, serum, or other additives may increase or decrease the aqueous solubility.

**Biological Activity:** SAHA inhibits HDAC1 with IC<sub>50</sub> of 40.6 nM using the BPS Bioscience Fluorogenic HDAC Assay Kit (Cat. # 50033).

### HDAC 1



**Storage/Stability:** Store at or below -20°C.

**Quality Control:** The purity was determined by HPLC.

#### References:

1. Marks, P.A. *Oncogene* **26**: 1351 – 1356 (2007).
2. Cang, S., *et al.*, *J. Hematol. Oncol.* **2**: 22 (2009).
3. Zhang, C., *et al.*, *J. Invest. Dermatol.* **125**: 1045 – 1052 (2005).