

Suberohydroxamic Acid (SBHA)

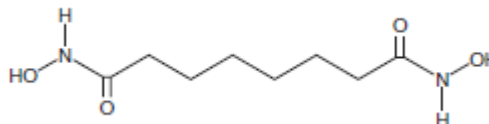
Catalog #: 27210

Lot #: 120529

Size: 100 mg

Structure:

CAS Registry #: 38937-66-5



Purity: ≥98%

Chemical Formula: C₈H₁₆N₂O₄

Molecular Weight: 204.2

Description: Suberohydroxamic acid (SBHA) is a competitive HDAC inhibitor. SBHA has been shown to cause cell differentiation, cell cycle arrest, and apoptosis.

Appearance: A crystalline solid

Solubility: Soluble in DMSO and DMF, purged with an inert gas. Solubility in DMSO and DMF is approximately 1.3 mg/ml and 0.1 mg/ml respectively. For maximum solubility in aqueous buffers, dissolve SBHA in DMSO then dilute with the aqueous buffer of choice. Do not store aqueous solutions for more than one day.

Biological Activity: SBHA inhibits HDAC1 with an IC₅₀ = 0.25 μM and HDAC3 with an IC₅₀ = 0.3 μM.

Storage/Stability: Store at or below -20°C for up to two years.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Richon, VM, *et al.*, *Proc. Natl. Acad. Sci USA*. 1998; **95**: 3003-3007.
2. Brinkmann, H., *et al.*, *J. Biol. Chem.* 2001; **276(25)**: 22491-22499.
3. Ning, L., *et al.*, *Oncologist*. 2008; **13**: 98-104.
4. Zhang, XD, *et al.*, *Mol. Cancer Ther.* 2004; **3(4)**: 425-435.
5. Zhang, XD, *et al.*, *Biochem. Pharmacol.* 2003; **66**: 1537-1545.
6. Flis, S., *et al.*, *Eur. J. Pharmacol.* 2010; **627**: 26-32.
7. Flis, S., *et al.*, *Biochem. Biophys. Res. Commun.* 2009; **387**: 336-341.