

CBHA

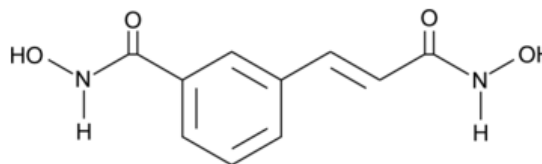
Catalog #: 27204

Lot #: 120208

Size: 5 mg

Structure:

CAS Registry #: 174664-65-4



Purity: ≥98%

Chemical Formula: C₁₀H₁₀N₂O₄

Molecular Weight: 222.2

Description: M-Carboxycinnamic Acid bishydroxamide, or CBHA, is a potent histone deacetylase (HDAC) inhibitor. CBHA has been found to induce apoptosis in nine different neuroblastoma cell lines in culture. It can also completely suppress neuroblastoma tumor growth in SCID mice at 200 mg/kg.

Appearance: A crystalline solid

Solubility: Soluble in DMSO and DMF at 20 mg/ml. First, dissolve in DMSO then dilute with the aqueous buffer of choice. Do not store in aqueous solution for more than one day.

Biological Activity: CBHA exhibits an IC₅₀ of 0.01 μM and 0.07 μM *in vitro* for HDAC1 and HDAC3 respectively.

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, V.M., *et al.*, *Proc. Natl. Acad. Sci.* 1998; **95**: 3003 – 3007.
2. Glick, R.D., *et al.*, *Cancer Res.* 1999; **59**: 4392 – 4399.
3. Coffey, D.C., *et al.*, *Cancer Res.* 2001; **61**: 3591 – 3594.