

HNHA

Catalog #: 27206

Lot #: 120403

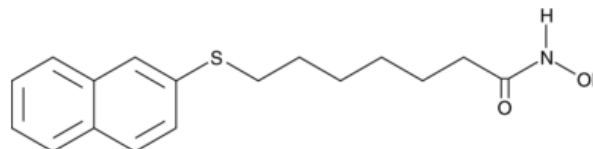
Size: 10 mg

Structure:

CAS Registry #: 926908-04-5

Purity: ≥98%

Chemical Formula: C₁₇H₂₁NO₂S



Molecular Weight: 303.4

Description: HNHA is a cell-permeable inhibitor of histone deacetylase (HDAC) activity. HNHA induces histone hyperacetylation and p21 transcription with inhibition of cell cycle progression in human fibrosarcoma HT1080 cells. In *in vivo* murine xenographs, HNHA is equally as effective as SAHA in inhibiting tumor growth. HNHA inhibits HUVEC growth and prevents tube formation and migration in response to VEGF. Also HNHA blocks retinal neovascularization and choroidal angiogenesis in mice.

Appearance: A crystalline solid

Solubility: Soluble in ethanol, DMSO and dimethyl formamide at 20 mg/ml. Sparingly soluble in aqueous buffers. First dissolve in DMSO, then dilute with aqueous buffer of choice. Do not store aqueous solutions for more than one day.

Biological Activity: HDAC is inhibited by HNHA with IC₅₀ = 100 nM.

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

Quality Control: The purity was determined by HPLC analysis.

References:

1. Kim, D.H., *et al.*, *Biochem. Biophys., Res. Commun.* 2007; **356**: 233 -328.
2. Kim, J.H., *et al.*, *Molecular Pharmaceutics* 2009; **6(2)**: 513-519.