

CI-994

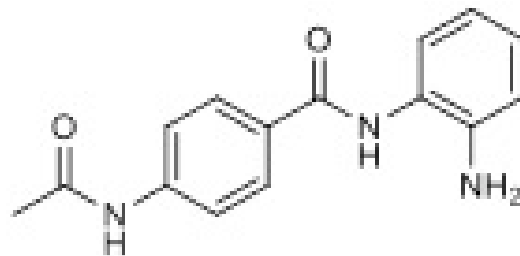
**Catalog No:** 14092, 14093**Format:** 25 mg, 5 mg**Chemical Properties:**

MW = 269.3

C<sub>15</sub>H<sub>15</sub>N<sub>3</sub>O<sub>2</sub>

CAS 112522-64-2

Physical Properties: Off-white powder

**Names:** N-acetyl-dinaline; 4-(Acetylamino)-N-(2-aminophenyl)benzamide; PD-123654

**Pharmacology:** Orally active histone deacetylase (HDAC) inhibitor (IC<sub>50</sub> = 0.9, 0.9, 1.2, and >20 μM for HDAC1, HDAC2, HDAC3 and HDAC8 respectively) (ref1). Mediates G<sub>1</sub> cell cycle arrest, inhibits proliferation and induces apoptosis *in vitro* and *in vivo* (ref 2,3). Increase neuroplasticity during memory extinction (ref 4). Protects beta cells from cytokine-induced apoptosis (ref 4).

**Solubilization:** May be dissolved in DMSO (25 mg/ml)**Fluorescent Properties:** N/A**Quality Control:**

&gt;98% (TLC); NMR (Conforms)

## References:

1. OM Moradei *et al. J. Med. Chem.* 2007, 50:5543
2. T Beckers *et al. Int. J. Cancer* 2007, 121:1138
3. M Loprevite *et al. Oncol. Res.* 2005, 15:39
4. DH Chou *et al. Chem. Biol.* 2012, 19:669

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.