

## Trichostatin A

**Catalog No:** 14038, 14039

**Format:** 5 mg, 1 mg

**Chemical Properties:**

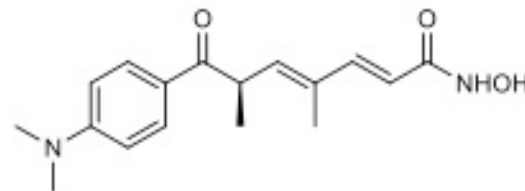
MW = 302.4

C<sub>17</sub>H<sub>22</sub>N<sub>2</sub>O<sub>3</sub>

CAS 58880-19-6

Physical Properties: Tan powder, MP = 150-151°C

**Names:** Trichostatin A, (2E,4E,6R)-7-(4-(Dimethylamino)phenyl)-N-hydroxy-4,6-dimethyl-7-oxo-2,4-heptadienamamide, TSA



**Pharmacology:** Potent and selective histone deacetylase (HDAC) inhibitor (K<sub>i</sub> = 3.4 nM)(ref 1). Induces reversion of ras-transformed cells to normal morphology (ref 2). Induces dedifferentiation of primordial germ cells into embryonic germ cells (ref 3). Cell permeable. Active *in vivo*.

**Solubilization:** May be dissolved in DMSO (15 mg/ml); Ethanol (3 mg/ml)

**Fluorescent Properties:** N/A

**Quality Control:**

>99% (HPLC); NMR (Conforms)

**References:**

1. M Yoshida *et al.* *J. Biol. Chem.* 1990, 265:17174
2. M Futamura *et al.* *Oncogene* 1995, 10:1119
3. G Durcova-Hills *et al.* *PLoS-One* 2008, 3:e3531

**Storage and Guarantee:** Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.