

## Panobinostat

**Catalog No:** 14044, 14045

**Format:** 25 mg, 5 mg

**Chemical Properties:**

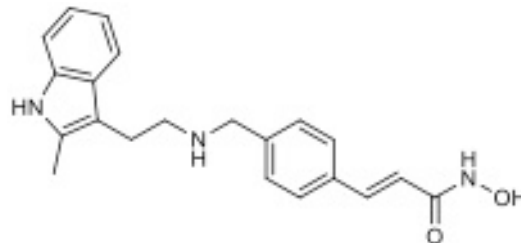
MW = 349.4

C<sub>21</sub>H<sub>23</sub>N<sub>3</sub>O<sub>2</sub>

CAS 404950-80-7

Physical Properties: Off white powder

**Names:** Panobinostat, (E)-N-hydroxy-3-(4-((2-(2-methyl-1H-indol-3-yl)ethylamino)methyl)phenyl)acrylamide



**Pharmacology:** A potent inhibitor of class I and II HDACs (ref 1). Cotreatment with panobinostat and an HSP90 inhibitor caused synergistic apoptosis in human CML-BC and AML cells (ref 2). A potent antimyeloma agent that overcomes drug resistance (ref 3).

**Solubilization:** May be dissolved in DMSO (100 mg/ml); Ethanol (5 mg/ml warm)

**Fluorescent Properties:** N/A

**Quality Control:**

≥98% (TLC); NMR (Conforms)

**References:**

1. L Geng *et al. Cancer Res.* 2006, 66:11298
2. P George *et al. Blood* 2005, 105:1768
3. P Maiso *et al. Cancer Res.* 2006, 66:5781

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