

Panobinostat

Catalog #: 27020

Size: 5 mg

CAS Registry #: 49843-98-3

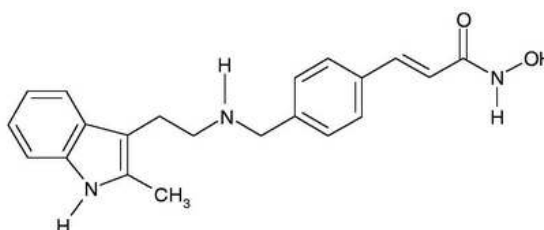
Purity: >99%

Chemical Formula: C₂₁H₂₃N₃O₂

Molecular Weight: 349.4

Lot #: 130925

Structure:



Description: Panobinostat is a histone deacetylase inhibitor. Inhibition of histone deacetylase (HDAC) by panobinostat is being researched in studies conducted for cutaneous T cell lymphoma, prostate cancer, myelodysplastic syndromes, breast cancer, chronic myelogenous leukemia, and other types of malignant disease.

Appearance: Yellow Solid.

Solubility: Soluble in DMSO at 200 mg/mL and ethanol 5 mg/mL. Poorly soluble in H₂O.

Biological Activity: Panobinostat was shown to inhibit several human multiple myeloma cell lines with an IC₅₀ of ≥ 20nM. Panobinostat was also shown to inhibit the proliferation of 7 of 8 tested pancreatic cell lines with a mean IC₅₀ of 0.09 μM.

Storage/Stability: Store at or below -20°C.

Quality Control: The purity was determined by HPLC analysis.

Reference(s):

1. Catley, L., *et al. Blood*. 2006 **108**: 3441-3449
2. Haefner, M., *et al. World J. Gastroenterol*. 2008 **14**:3681-3692