Tubastatin A hydrochloride



Catalog No: 14084, 14085

Format: 25 mg, 5 mg

Chemical Properties:

MW = 371.9

C₂₀H₂₁N₃O₂• HCI

CAS 1310693-92-5

Physical Properties: White powder

Names: Tubastatin A, N-Hydroxy-4-((2-methyl-3,4-dihydro-1H-pyrido[4,3-b]-

indol-5(2H)-yl)methyl)benzamide, hydrochloride

Pharmacology: Potent and selective histone deacetylase 6 (HDAC 6) inhibitor (IC $_{50}$ = 15 nm; >1000-fold selective against other HDACs and 60-fold for HDAC8))(ref 1). Inhibits TNF α (IC $_{50}$ = 272 nm) and IL-6 (IC $_{50}$ = 712 nm) in LPS-stimulated human macrophages and displays anti-inflammatory activity in animal models (ref 2). Improves survival in animal models of sepsis (ref 3). Improves cognition in Alzheimers disease transgenic mice (ref 4). Cell permeable.

Solubilization: May be dissolved in DMSO (50 mg/ml)

Fluorescent Properties: N/A

Quality Control:

99% (HPLC); NMR (Conforms)

References:

1. KV Butler et al. J. Am. Chem. Soc. 2010, 132:10842

2. S Vishwakarma et al. Int. Immunopharmacol. 2013, 16:72

3. T Zhao et al. Surg. Res. 2014, 190:647

4. L Zhang et al. J. Alzheimers Dis. 2014, 41:1193

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.

HCI NHOH

Chemical structure of Tubastatin A hydrochloride.