Romidepsin



Catalog No: 14082, 14083

Format: 5 mg, 1 mg

Chemical Properties:

MW = 540.7 $C_{24}H_{36}N_4O_6S_2$ CAS 128517-07-7

Physical Properties: Colorless lypohilized wax (vials appear empty) **Names:** Romidepsin, FK228, FK 228, FR901228, Depsipeptide

Pharmacology: A potent and selective inhibitor of class I histone deacetylases (HDACs), IC_{50} = 36, 47, 510 and 14,000 nM for HDAC1, 2, 4 and 6 respectively (ref 1). Induces apoptosis in a variety of cell lines (ref 2) and displays anti-tumor activity in mouse models (ref 3). Recently approved for clinical uses in T-cell lymphoma (ref 4).

Solubilization: May be dissolved in DMSO (10 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (HPLC); NMR (Conforms)

References:

1. R Furumai et al. Cancer Res. 2002, 62:4916

2. J Panicker et al. Cell Cycle 2010, 9:1830

3. H Ueda et al. J. Antibiot. (Tokyo) 1994, 47:315

4. MK VanderMolin et al. J. Antibiot. (Tokyo) 2011, 64:525

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 for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.

