

Romidepsin

Catalog No: 14083

Format: 1 mg

Chemical Properties:

MW = 540.7

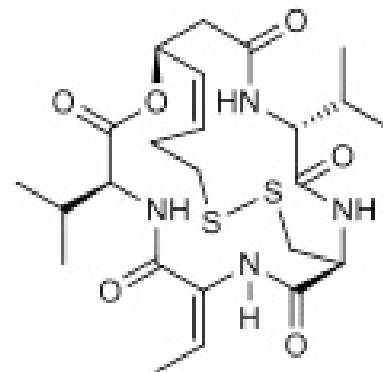
C₂₄H₃₆N₄O₆S₂

CAS 128517-07-7

Physical Properties: Colorless lyophilized wax (vials appear empty)

Names: Romidepsin; FK228; FR901228; Depsipeptide

Pharmacology: A potent and selective inhibitor of class I histone deacetylases (HDACs), IC₅₀ = 36, 47, 510 and 14,000 nM for HDAC1, 2, 4 and HDAC6 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).



Chemical structure of Romidepsin.

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 guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.