

Mocetinostat

Catalog No: 14136, 14137

Format: 25 mg, 5 mg

Chemical Properties:

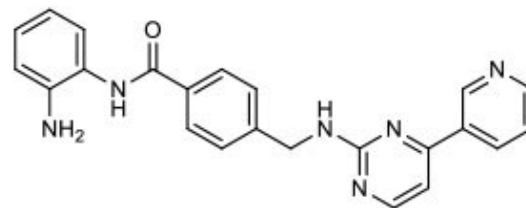
MW = 396.4

C₂₃H₂₀N₆O

CAS 726169-73-9

Physical Properties: Off-white crystalline powder

Names: Mocetinostat; N-(2-aminophenyl)-4-[(4-pyridin-3-ylpyrimidin-2-ylamino)methyl]benzamide; MGCD0103



Chemical structure of Mocetinostat.

Pharmacology: Class I, isoform-selective HDAC inhibitor (IC₅₀ = 0.15, 0.29, 1.66 and 0.59 μM for HDAC1, 2, 3 and 11 respectively) (ref 1). Induces hyperacetylation of histones, induces expression of the tumor suppressor p21WAF1 and inhibits proliferation of human cancer cells (ref 2). Displays antifibrotic effects in ischemic heart failure (ref 3). Attenuates the development of hypersensitivity in models of neuropathic pain (ref 4). Active *in vivo* (ref 5).

Solubilization: May be dissolved in DMSO (25 mg/ml)

Fluorescent Properties: N/A

Quality Control:

>98% (HPLC); NMR (Conforms)

References:

1. N Zhou *et al. J. Med. Chem.* 2008, 51:4072
2. S Raeppl *et al. Bioorg. Med. Chem. Lett.* 2009, 19:644
3. H Nural-Govener *et al. Int. J. Mol. Sci.* 2015, 16:11482
4. F Denk *et al. Pain* 2013, 154:1668
5. C Bonfils *et al. Clin. Cancer Res.* 2008, 14:3441

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 2 years. Store solutions at -20°C for up to 2 months. 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.