CUDC-101



Catalog No: 14060, 14061

Format: 25 mg, 5 mg

Chemical Properties:

MW = 434.5

 $C_{24}H_{26}N_4O_4$

CAS 1012054-59-9

Physical Properties: White to off-white powder; MP = 188-190°C

Names: CUDC101, 7-((4-((3-ethynylphenyl)amino)-7-methoxyquinazolin-6-yl)

oxy)-N-hydroxyheptanamide

Pharmacology: A novel hybrid dual-acting HDAC and receptor tyrosine kinase inhibitor. It is a potent HDAC inhibitor which also inhibits EGFR and HER2 (IC $_{50}$ = 4.4 nM, 2.4 nM and 15.7 nM respectively) (ref 1). It not only blocks EGFR and HER2 but also attenuates multiple compensatory pathways such as AKT, HER3 and MET which enable tumor cells to escape the effects of conventional EGFR/HER2 inhibitors (ref 1,2).

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

Fluorescent Properties: N/A

Quality Control:

>99% (HPLC); NMR (Conforms)

References:

- 1. CJ Lai et al. Cancer Res. 2010, 70:3647
- 2. X Cai et al. J. Med. Chem. 2010, 53:2000

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 arrival.

