

## CUDC-101

Epigenetic Enzyme Inhibitor

### Catalog # H83-904

Lot # P3320-13

CAS # 1012054-59-9

### Product Description

Molecular Formula: C<sub>24</sub>H<sub>26</sub>N<sub>4</sub>O<sub>4</sub>

Appearance: White to off-white powder

Molecular Weight: 434.5

Purity: >99% (HPLC); NMR (Conforms)

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

### Alias

7-((4-((3-ethynylphenyl)amino)-7-methoxyquinazolin-6-yl)oxy)-N-hydroxyheptanamide

### Specific Activity

Dual HDAC & EGFR Inhibitor

### Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

Store solutions at -20°C for up to 1 month.

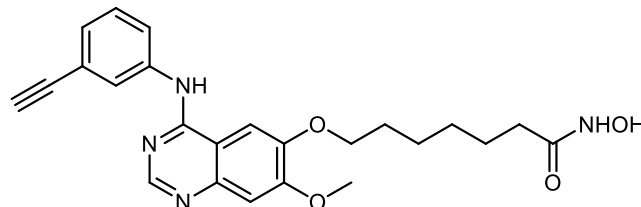
### Scientific Background

A novel hybrid dual-acting HDAC and receptor tyrosine kinase inhibitor. It is a potent HDAC inhibitor which also inhibits EGFR and HER2, IC<sub>50</sub> = 4.4, 2.4 and 15.7 nM respectively<sup>1</sup>. 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<sup>1,2</sup>.

### References

1. C.J. Lai et al. Cancer Res. 2010 70:3647
2. X. Cai et al. J. Med. Chem. 2010 53:20

### Molecular Structure



## CUDC-101

Epigenetic Enzyme Inhibitor

Catalog #	H83-904
CAS #	1012054-59-9
Lot #	P3320-13
Purity	>99% (HPLC)
Format	White to off-white powder
Solubilization	May be dissolved in DMSO (25 mg/ml)
Storage & Shipping	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**