

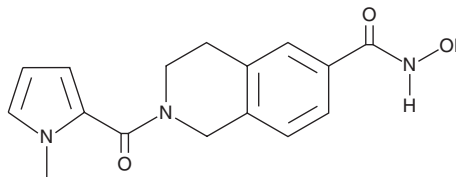
# PRODUCT INFORMATION



## HDAC6 Inhibitor

Item No. 15200

**CAS Registry No.:** 1259296-46-2  
**Formal Name:** 1,2,3,4-tetrahydro-N-hydroxy-2-[(1-methyl-1H-pyrrol-2-yl)carbonyl]-6-isoquinolinecarboxamide  
**Synonym:** Histone Deacetylase 6  
**MF:** C<sub>16</sub>H<sub>17</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 299.3  
**Purity:** ≥95%  
**Supplied as:** A solution in acetonitrile  
**Storage:** -20°C  
**Stability:** ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

HDAC6 inhibitor is supplied as a solution in acetonitrile. To change the solvent, simply evaporate the acetonitrile under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of HDAC6 inhibitor in these solvents is approximately 30 mg/ml.

### Description

HDAC6 is a predominantly cytoplasmic enzyme that targets  $\alpha$ -tubulin, cortactin, and heat shock protein 90, as well as other substrates.<sup>1</sup> In this way, it impacts development, proliferation, invasion, and tumorigenesis.<sup>1</sup> HDAC6 Inhibitor is a potent and selective inhibitor of HDAC6 (IC<sub>50</sub> = 36 nM) that poorly blocks other HDAC enzymes.<sup>2</sup> It is cell permeable, inhibiting the acetylation of tubulin in cells with an IC<sub>50</sub> value of 210 nM.<sup>2</sup>

### References

1. Yang, P.H., Zhang, L., Zhang, Y.J., *et al.* HDAC6: Physiological function and its selective inhibitors for cancer treatment. *Drug Discov. Ther.* **7(6)**, 233-242 (2013).
2. Blackburn, C., Barrett, C., Chin, J., *et al.* Potent histone deacetylase inhibitors derived from 4-(aminomethyl)-N-hydroxybenzamide with high selectivity for the HDAC6 isoform. *J. Med. Chem.* **56(18)**, 7201-7211 (2013).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 08/16/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)