

# PCI-34051

## HDAC8 inhibitor

PCI-34051 is a potent and selective HDAC8 inhibitor with greater than 200-fold selectivity over HDACs 1, 2, 3, 6, 10. Unlike broad-spectrum HDAC inhibitors, it does not cause detectable histone or tubulin acetylation. It is a very useful tool for exploring the role of HDAC8 in cellular function.

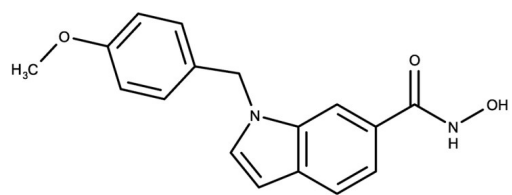
### Ordering Information

[Order Online »](#)

ENZ-CHM445-0025	25mg
-----------------	------

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** N-Hydroxy-1-[(4-methoxyphenyl)methyl]-1H-indole-6-carboxamide

**Appearance** Off-white solid.

**CAS** 950762-95-5

**Couple Type** Inhibitor

**Formula**  $C_{17}H_{16}N_2O_3$

**Identity** Determined by NMR.

**MW** 296.33

**Purity** ≥98% (TLC)

**Solubility** Soluble in DMSO (50 mg/ml).

**Last modified: January 24, 2025**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)