

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.

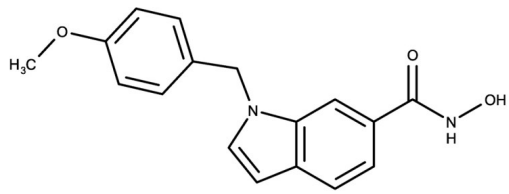
Ordering Information

[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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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