

MC 1293

HDAC1 inhibitor

Inhibitor of histone deacetylase 1 (HDAC1) ($IC_{50} = 4.5\mu M$) and maize histone deacetylase HD2 ($IC_{50} = 1.9\mu M$).

Citations: 3

[View Online »](#)

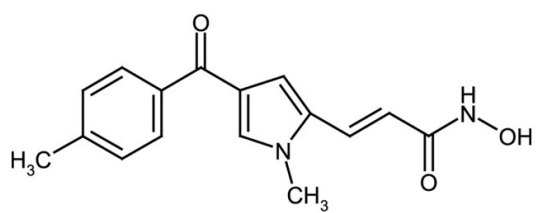
Ordering Information

[Order Online »](#)

ALX-270-344-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-(4-Toluoyl-1-methyl-1H-2-pyrrolyl)-N-hydroxy-2-propenamide
Appearance	Off-white powder.
Couple Target	HDAC, HDAC1, HDAC2
Couple Type	Inhibitor
Formula	$C_{16}H_{16}N_2O_3$
Identity	Determined by 1H-NMR.
MW	284.3
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

Last modified: May 29, 2024



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)