

M344

HDAC inhibitor

Potent inhibitor of histone deacetylases (HDACs) ($IC_{50} \leq 1 \mu M$).

Citations: 2

[View Online »](#)

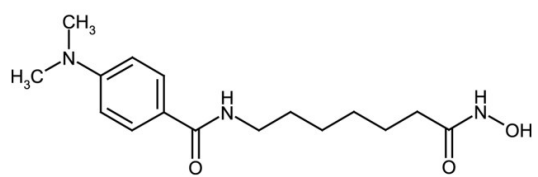
Ordering Information

[Order Online »](#)

ALX-270-297-M001	1mg
ALX-270-297-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	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	4-Dimethylamino-N-(6-hydroxycarbamoylhexyl)-benzamide
Appearance	White crystalline solid.
CAS	251456-60-7
Couple Target	HDAC
Couple Type	Inhibitor
Formula	$C_{16}H_{25}N_3O_3$
Identity	Determined by NMR.
MW	307.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (1mg/ml)

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)