

# M344

## HDAC inhibitor

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

Citations: 2

[View Online »](#)

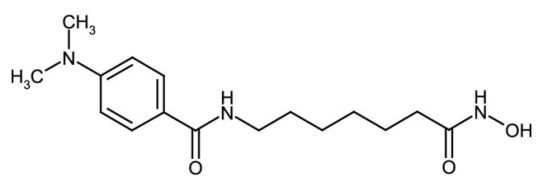
### Ordering Information

[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)