

# BML-210

## HDAC inhibitor

HDAC inhibitor developed at Enzo Life Sciences. Inhibits HDAC activity in HeLa nuclear extracts ( $IC_{50}$ =5-10 $\mu$ M).

Citations: 3

[View Online »](#)

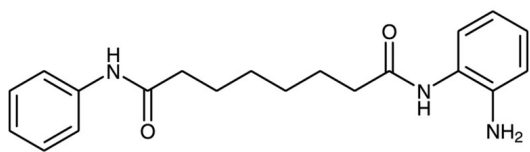
## Ordering Information

[Order Online »](#)

BML-GR330-0001	1mg
----------------	-----

## 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

**Appearance** Tan solid.

**CAS** 537034-17-6

**Couple Target** HDAC

**Couple Type** Inhibitor

**Formula**  $C_{20}H_{25}N_3O_2$

**MW** 339.4

**Purity** ≥98% (TLC)

**Solubility** Soluble in DMSO (25mg/ml) or 100% ethanol (10mg/ml, warm)

**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-be@enzolifesciences.com](mailto:info-be@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)