

Apicidin

HDAC inhibitor

Potent inhibitor of histone deacetylase (HDAC). Inhibits proliferation. Induces cell cycle arrest at the G1 phase. At 100nM it induces a long lasting hyperacetylation of histone H4 while that induced by trichostatin is transient. Stimulates apoptosis. Apoptosis is induced via induction of Fas/Fas ligand. Displays potent antiangiogenic effects and dramatically decreases HIF-1 α protein levels and transcriptional activity in human and mouse tumor cell lines. Antiprotozoal.

Citations: 20

[View Online »](#)

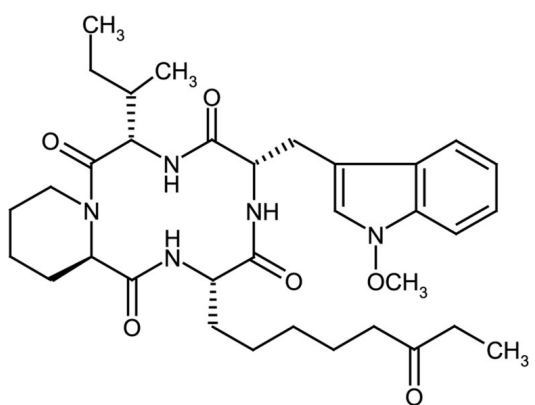
Ordering Information

[Order Online »](#)

BML-GR340-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	cyclo-L-(2-Amino-8-oxodecanoyl)-L-(N-methoxy-tryptophan)-L-isoleucyl-D-pipecolinyl, Cyclo(N-O-methyl-L-tryptophanyl-L-isoleuciny-L-D-pipecolinyl-L-2-amino-8-oxodecanoyl)
Appearance	White powder.
CAS	183506-66-3
Couple Target	HDAC
Couple Type	Inhibitor
Formula	$C_{34}H_{49}N_5O_6$
MW	623.8
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO or 100% ethanol.



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)