

Actinonin

Protease inhibitor

Antibiotic. Inhibitor of aminopeptidase M, N and leucine and alanyl aminopeptidase and peptide deformylase (PDF). Inhibits MMP-1 (K_i =300nM), MMP-2 (IC_{50} ~90nM at MMP-2=5nM), MMP-3 (K_i =1'700nM), MMP-7 (IC_{50} ~3'500nM at MMP-7=19nM), MMP-8 (K_i =160nM), MMP-9 (K_i =330nM), MMP-10 (IC_{50} ~500nM at MMP-10=22nM), MMP-12 (IC_{50} ~350nM at MMP-12=4nM), and MMP-13 (IC_{50} ~100nM at MMP-13=13nM), but not MMP-14 (IC_{50} >1mM at MMP-14=13.6nM). Shows antitumor activity *in vitro* and *in vivo*. Apoptosis inducer.

Citations: 14

[View Online »](#)

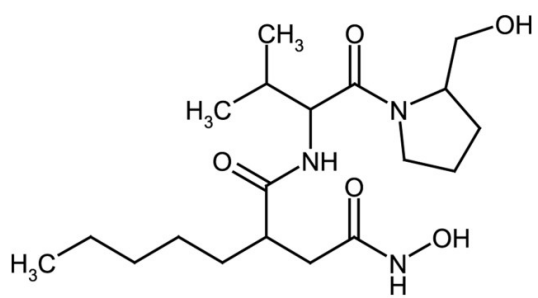
Ordering Information

[Order Online »](#)

ALX-260-128-M005	5mg
------------------	-----

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

Alternative Name	3[[1-[[2-(Hydroxymethyl)-1-pyrrolidinyl]-carbonyl]-2-methylpropyl]-carbamoyl]-octanohydroxamic acid
Appearance	White to off-white powder.
CAS	13434-13-4
Couple Target	Aminopeptidase, MMP, MMP-1, MMP-10, MMP-12, MMP-13, MMP-2, MMP-3, MMP-7, MMP-8, MMP-9, Peptide deformylase
Couple Type	Inhibitor
Formula	$C_{19}H_{35}N_3O_5$
MW	385.5
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol (20mg/ml) or DMSO (>20mg/ml).
Source	Synthetic.

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

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)