

Actinonin

Protease inhibitor

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

Citations: 14

[View Online »](#)

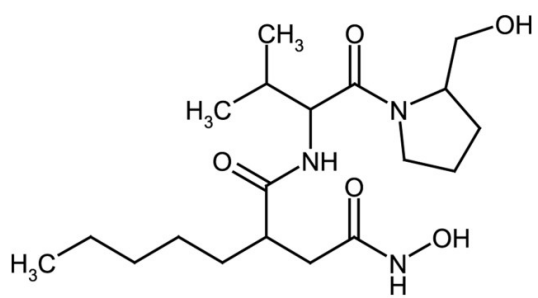
Ordering Information

[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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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