

Pepstatin A (synthetic)

Protease inhibitor

Inhibitor of aspartate (acid) proteases, including pepsin, cathepsin D, renin, chymosin, bacterial aspartic proteinases and HIV proteases. Efficiently inhibits Fcγ receptor induction by IFN-γ. Blocks the endocytic pathway followed by an antibody when it enters the cell at the level of early endosomal compartments, resulting in inhibition of autoantigen presentation or viral transmission. Suppresses p-53-dependent apoptosis in lymphoid cells as well as TNFα-induced apoptosis in U937 cells. Inhibits degradation of autophagic cargo inside autophagolysosomes.

Citations: 25

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-085-M005	5mg
ALX-260-085-M025	25mg
ALX-260-085-M100	100mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Iva-Val-Val-Sta-Ala-Sta, i-Valeryl-L-Val-L-Val-AHMHA-L-Ala-AHMHA
Appearance	White to off-white solid.
CAS	26305-03-3
Formula	$C_{34}H_{63}N_5O_9$
MI	14: 7147
MW	685.9
Purity	≥98% (HPLC)
RTECS	SC6155000
Sequence	i-Valeryl-L-Val-L-Val-AHMHA-L-Ala-AHMHA (AHMHA = (3S,4S)-4-Amino-3-hydroxy-6-methyl-heptanoic acid)
Solubility	Soluble in DMSO (10mg/ml), methanol (1mg/ml) or acetic acid.

Source

Synthetic.



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-
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)

Last modified: May 29, 2024