

# Pepstatin A (synthetic)

## Protease inhibitor

Inhibitor of aspartate (acid) proteases, including pepsin, cathepsin D, renin, chymosin, bacterial aspartic proteinases and HIV proteases. Efficiently inhibits Fc $\gamma$  receptor induction by IFN- $\gamma$ . 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 TNF $\alpha$ -induced apoptosis in U937 cells. Inhibits degradation of autophagic cargo inside autophagolysosomes.

Citations: 25

[View Online »](#)

## Ordering Information

[Order Online »](#)

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

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

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

Last modified: May 29, 2024