

# 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<br>(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-  
eu@enzolifesciences.com](mailto:info-eu@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)