

# Amastatin . hydrochloride

## Aminopeptidase inhibitor

Competitive inhibitor of human serum aminopeptidase A, pig kidney leucine aminopeptidase, leucine aminopeptidase, tyrosine aminopeptidase, but not of aminopeptidase B. Has been used to potentiate the action of some bioactive peptides by inhibiting their degradation by aminopeptidases.

Citations: 5

[View Online »](#)

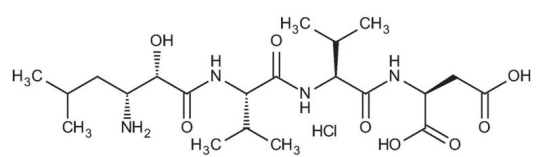
## Ordering Information

[Order Online »](#)

ALX-260-003-M001	1mg
ALX-260-003-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** [(2S,3R)-3-Amino-2-hydroxy-5-methyl-hexanoyl]-Val-Val-Asp-OH . HCl

**Appearance** White to off-white powder.

**CAS** 100938-10-1

**Formula**  $C_{21}H_{38}N_4O_8 \cdot HCl$

**MW** 474.6 . 36.5

**Purity** ≥98% (HPLC)

**Solubility** Soluble in 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)