# 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 »

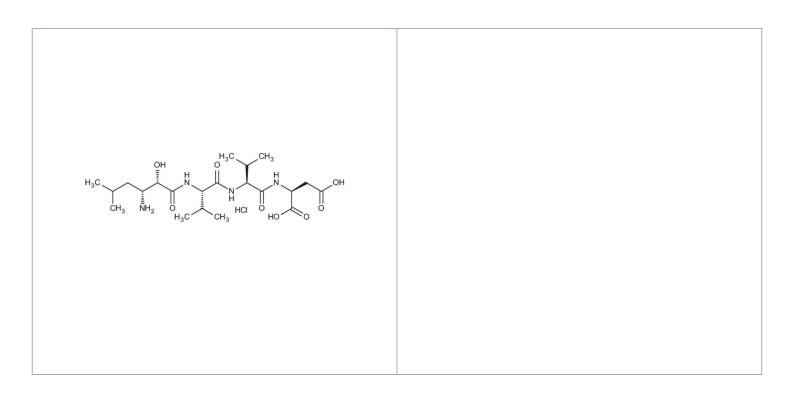
## **Ordering Information**

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 . HCI

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

**CAS** 100938-10-1

**Formula**  $C_{21}H_{38}N_4O_8$ . HCI

474.6 . 36.5 MW

**Purity** ≥98% (HPLC)

**Solubility** Soluble in acetic acid.

Synthetic. Source

Last modified: May 29, 2024

uk@enzolifesciences.com