

# Antipain

## Papain inhibitor

Inhibits papain and trypsin and to a small extent plasmin. Antipain is more specific for papain and trypsin than is leupeptin (Prod. No. ALX-260-009). The inhibitory potency of antipain is 100-fold higher than that of elastatinal. Acts by forming a hemiacetal adduct between its aldehyde group and the active serine of the proteinase.

Citations: 3

[View Online »](#)

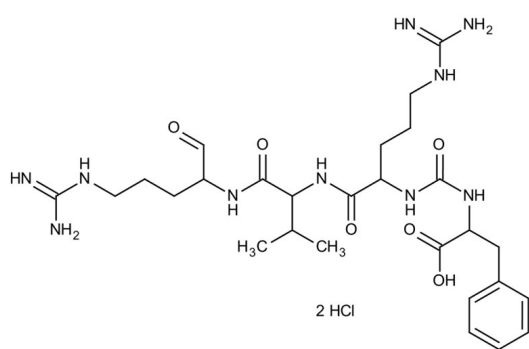
## Ordering Information

[Order Online »](#)

ALX-260-093-M005	5mg
ALX-260-093-M025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

**Use/Stability** As indicated on product label or CoA when stored as recommended.

**Handling** Protect from light.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Activity** Inhibitory Activity  $\geq 30$  U/mg (Trypsin, L-BAPA)

**Alternative Name** [(S)-1-Carboxy-2-phenylethyl]-carbamoyl-Arg-Val-arginal . 2HCl

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

**CAS** 37682-72-7

**Couple Target** Papain

**Couple Type** Inhibitor

**Formula**  $C_{27}H_{44}N_{10}O_6 \cdot 2HCl$

**MW** 604.7 . 72.9

**Purity**  $\geq 97\%$  (HPLC)

**Sequence** [(S)-1-Carboxy-2-phenylethyl]-carbamoyl-Arg-Val-arginal

**Solubility** Soluble in water, methanol, 100 % ethanol or acetonitrile.

**Source** Synthetic.