

Chymostatin

Protease inhibitor

Inhibitor of serine proteinases having a chymotrypsin-like specificity like α -, β -, δ - and γ -chymotrypsin, chymases, cathepsin G, and most cysteine proteinases including papain and cathepsin A, B, H, and L. Mixture of type A (R=L-Leu), type B (R=L-Val), type C (R=L-Ile).

Citations: 1

[View Online »](#)

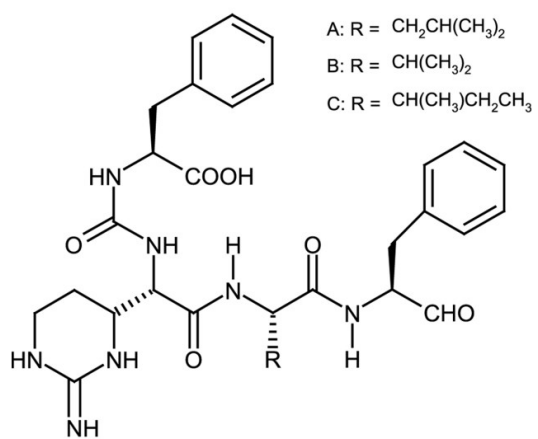
Ordering Information

[Order Online »](#)

ALX-260-005-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 (S)-1-Carboxy-2-phenylethyl)-carbamoyl- α -[2-iminohehexahydro-4(S)-pyrimidyl]-(S)-glycyl-X-phenylalaninal

Appearance White to off-white powder.

CAS 9076-44-2

Couple Target Cathepsin, Chymase, Chymotrypsin, Papain

Couple Type Inhibitor

Formula $C_{31}H_{41}N_7O_6$

MW 607.7

Purity $\geq 95\%$ TLC (mixture of A, B, C)

Solubility Soluble in DMSO (20mg/ml) or glacial acetic acid.

Source Microbial.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-be@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com