

Apstatin

Aminopeptidase inhibitor

Apstatin is a potent and selective inhibitor of aminopeptidase P (APP), $K_i = 2.6 \mu\text{M}$ for purified rat lung membrane-bound APP. It blocks the APP-mediated degradation of bradykinin. Limits myocardial infarct size alone or with ACE inhibitors.

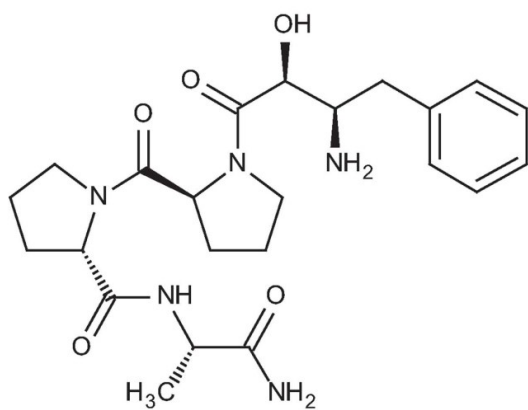
Ordering Information

[Order Online »](#)

BML-PI145-0001	1mg
----------------	-----

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-[(2S,3R)-3-Amino-2-hydroxy-4-phenylbutanoyl]-L-prolyl-L-prolyl-L-alaninamide

Appearance Solid.

CAS 160470-73-5

Formula $C_{23}H_{33}N_5O_5$

Formulation Lyophilized.

MW 459.5

Purity ≥95% (HPLC)

Sequence N-[(2S,3R)-3-Amino-2-hydroxy-4-phenylbutanoyl]-L-prolyl-L-prolyl-L-alaninamide

Solubility Soluble in water (>25 mg/ml) or DMSO (20 mg/ml).

Last modified: May 29, 2024



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)