

Z-prolyl-prolinal

Endopeptidase inhibitor

A potent and selective inhibitor of prolyl endopeptidase (PEP/PREP/POP, EC 3.4.21.26). $K_i=0.35$ nM for mouse brain enzyme. Inhibits the cleavage of arginine-vasopressin, gonadotropin-releasing hormone, and other peptides.

Citations: 4

[View Online »](#)

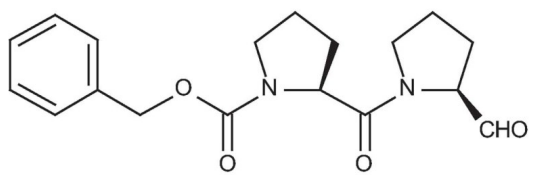
Ordering Information

[Order Online »](#)

BML-PI112-0005	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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Z-PP-CHO
Appearance	White solid.
CAS	86925-97-5
Formula	$C_{18}H_{22}N_2O_4$
MW	330.4
Purity	≥97% (HPLC)
Sequence	Z-Pro-Pro-CHO
Solubility	Soluble in DMSO.

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-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@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)