

POP/PREP (human), (recombinant)

POP/PREP is a cytoplasmic, membrane-associated, or serum serine protease that cleaves peptide hormones and neuropeptides such as substance P, oxytocin, neurotensin, angiotensins, and bradykinin. This enzyme is related to DPPIV and is useful for specificity screening of DPPIV inhibitors.

Ordering Information

[Order Online »](#)

BML-SE545-0010	10µg
-----------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Shipping

Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name

Prolyl oligopeptidase, Prolyl endopeptidase, Post-proline cleaving enzyme

Application Notes

Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.

MW

81.6 kDa

Source

Produced in *E. coli*.

UniProt ID

P48147

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)