

DPPIII (human), (recombinant)

DPPIII is a metalloproteinase that cleaves biologically active peptides such as enkephalins, angiotensins, and endomorphins, and has been implicated in cancer, pain, and inflammation.

Ordering Information

[Order Online »](#)

BML-SE529-0010	10µg
-----------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dipeptidyl Peptidase III, DPP3

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

MW 109 kDa

Source Produced in insect cells. Active recombinant human DPPIII with an N-terminal purification tag. Produced in a baculovirus expression system.

UniProt ID Q9NY33



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)