

DPP8 (human), (recombinant) (GST- tag)

The biological function of DPP8 is not known, but it is found in the cytoplasm and may have roles in cell adhesion, migration, and apoptosis. This enzyme is related to DPPIV and is useful for specificity screening of DPPIV inhibitors to avoid toxic side effects.

Ordering Information

[Order Online »](#)

BML-SE527-0010	10µg
-----------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	The enzyme is stable on ice for at least several hours. However, it is recommended that thawing and dilution of the enzyme be done within as short a time as possible before start of the assay. NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing could lead to loss of activity.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Dipeptidyl peptidase 8
Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
Formulation	Liquid. In 40mM TRIS-HCl, pH 8.0, containing 110mM sodium chloride, 2.2mM potassium chloride, and 20% glycerol.
GenBank ID	NM_197960
MW	129 kDa
Purity	≥70% (SDS-PAGE)
Source	Produced in insect cells. Active recombinant human DPP8 is fused at the N-terminus to a GST-tag. Produced in a baculovirus expression system.
UniProt ID	Q6V1X1



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-
fr@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)