

DPPIV (human), (recombinant)

Highly active

DPPIV is involved in diseases such as diabetes (*via* cleavage of GIP and GLP-1) and cancer.

Citations: 5

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE434-0025	25MU
----------------	------

Manuals, SDS & CofA

[View Online »](#)

kDa

148

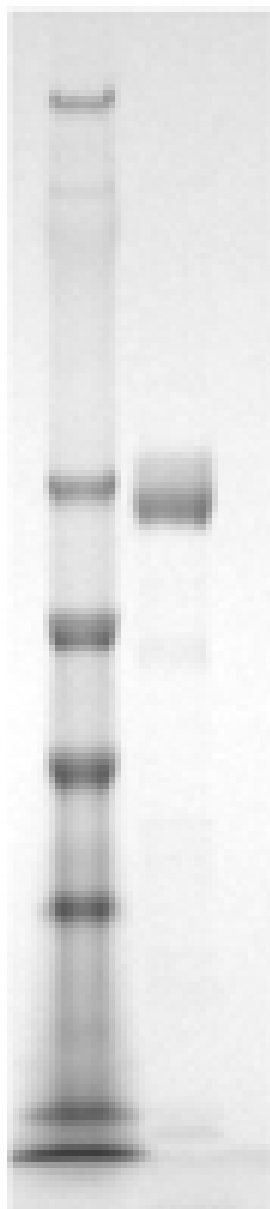
98

64

50

36

22



1

2

SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 1.0 μ g of purified Human DPPIV protein.

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. After initial defrost, the remaining, undiluted and unused enzyme should be refrozen quickly by, for example, snap-freezing in a dry ice ethanol bath or liquid nitrogen. Aliquot product shall be stored -80°C. Avoid repeated freeze/defrost cycles. 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.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Dipeptidyl Peptidase IV, DPP4, CD26
Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
Formulation	Liquid. In 10 mM TRIS, pH 8.0, 100mM NaCl.
MW	85.4 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in Sf9 insect cells using baculovirus expression system. Active recombinant human DPPIV (aa 29-766). This represents a naturally-occurring cleaved (soluble) form of DPPIV.
Specific Activity	One unit produces 1µmole free pNA per minute, using as substrate 100 µM H-Gly-Pro-pNA (Prod. No. BML-P188) in 50mM TRIS pH 7.5 at 37°C.
UniProt ID	P27487



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com