## 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** 

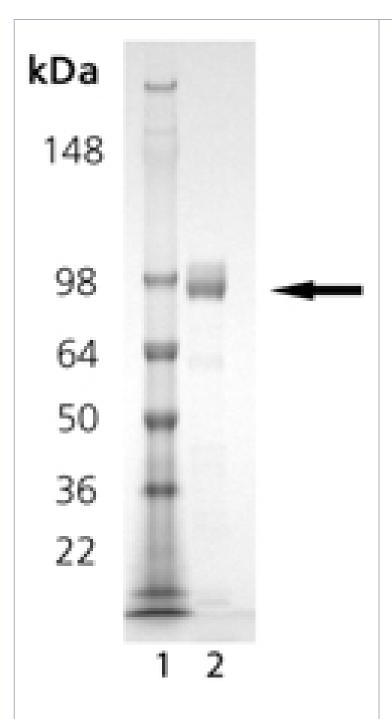
Order Online »

BML-SE434-0025

25MU

Manuals, SDS & CofA

**View Online** »



SDS-PAGE Analysis: Lane 1: MWM; Lane 2: 1.0  $\mu g$  of purifed 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com