DPP8 (human), (recombinant) (GSTtag)

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

