DPP9 (human), (recombinant)

The biological function of DPP9 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-SE528-0010 10

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 9

Application NotesUseful tool to study enzyme kinetics, cleave target

substrates, and screen for inhibitors.

Formulation Liquid. In 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM

KCI, 20% glycerol, and 3 mM DTT.

MW 124 kDa

Source Produced in insect cells. Active recombinant human DPP9

with an N-terminal purification tag. Produced in a

baculovirus expression system.

UniProt ID Q86TI2

Last modified: May 29, 2024

