

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.

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 Notes Useful 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 KCl, 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



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