

# 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	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 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



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

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

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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