

Data Sheet

DPP9

Human, recombinant, N-terminal GST tag

Catalog #: 80090

Lot#: 2000

Conc.: 2 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0,
124 mM NaCl, 2.4 mM KCl, 18 mM
Glutathione, 10% glycerol, and 3 mM DTT.

Stability: >6 months at -80°C

References:

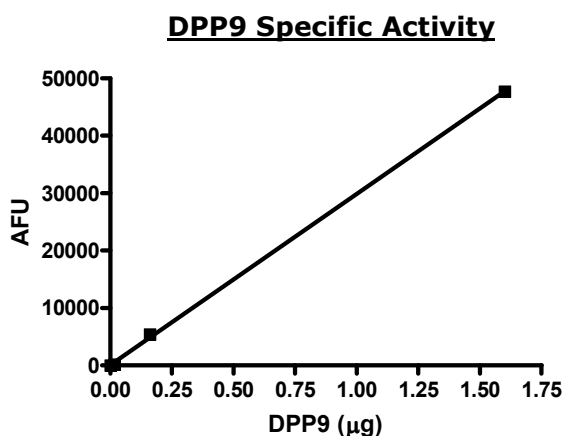
1. Bjelke, J.R. et al., Biochem. J. 396 (2), 391-399 (2006)
2. Yu, D.M. et al., FEBS J. 273 (11), 2447-2460 (2006)

Description: Human DPP9 (GenBank Accession No. NM_139159), (30-end) with N-terminal GST tag, MW=124 kDa, expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 536 U/ug. One U = 1 pmol/min, Assay condition: 25 mM Tris/Cl, pH7.5, 0.1% BSA, 5 uM substrate (H-Gly-Pro-AMC), and 3 ng/ul DPP9. Incubation condition: 1 min at room temperature.

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

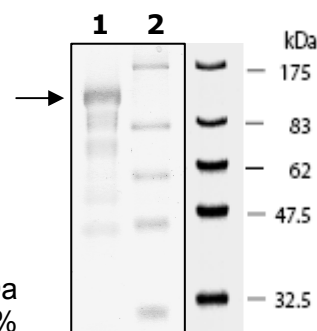
Quality Assurance



10% SDS-PAGE Coomassie staining

Lane 1:
7 µg DPP9
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 122.6 kDa
Purity: ≥70%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com