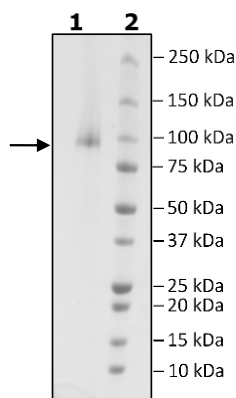


## Product Information

<b>Construct:</b>	DPP4 (29-766(end)-His)
<b>Concentration:</b>	0.63 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM Imidazole, 10% glycerol
<b>Expression System:</b>	Sf9
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_001935
<b>MW:</b>	87 kDa
<b>Purity:</b>	≥90%
<b>Specific Activity:</b>	≥4,000 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to the DPP4 Assay Kit (BPS Bioscience #80204) with various amounts of DPP4 enzyme.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

