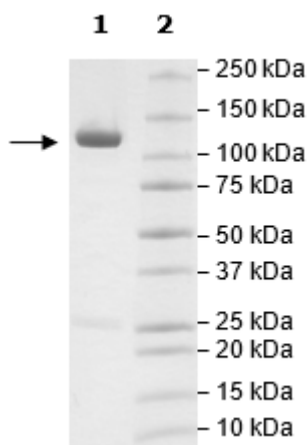


Product Information

Construct:	DPP9 (GST-1-863(end))
Concentration:	1.35 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT, and 16 mM glutathione
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001384623
MW:	124 kDa
Purity:	≥90%
Specific Activity:	250 pmol/min/μg
Assay Conditions:	Assay was done according to DPP9 Assay Kit (BPS Bioscience #80209) with DPP9 enzyme titrated at various concentrations and incubated at room temperature for 10 minutes.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



DPP9 Activity

