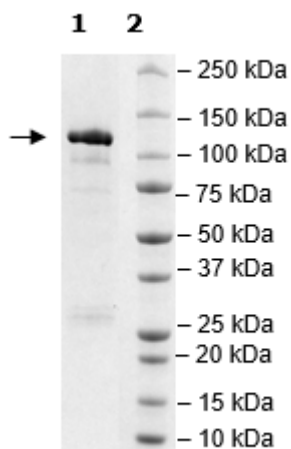


## Product Information

<b>Construct:</b>	DPP8 (GST-His-1-898(end))
<b>Concentration:</b>	2.96 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT
<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_197960
<b>MW:</b>	129 kDa
<b>Purity:</b>	80%
<b>Specific Activity:</b>	≥560 pmol/min/μg
<b>Assay Conditions:</b>	Assay was done according to the DPP8 Assay Kit (BPS Bioscience #80208) with DPP8 enzyme titrated from 80 ng/rxn-0 ng/rxn and incubated at room temperature for 10 minutes.
<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

