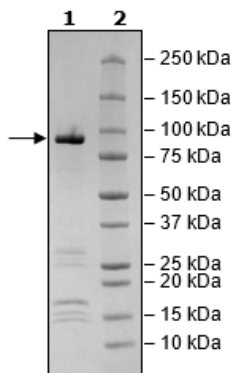


Product Information

Construct:	PARP2 (GST-2-583(end))
Concentration:	0.37 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, 5 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_005484
MW:	96 kDa
Purity:	72%
Assay Conditions:	Assay was performed according to PARP2 Chemiluminescent Assay Kit (BPS Bioscience #80552) with PARP2 serially diluted at various concentrations.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



PARP2 Activity

