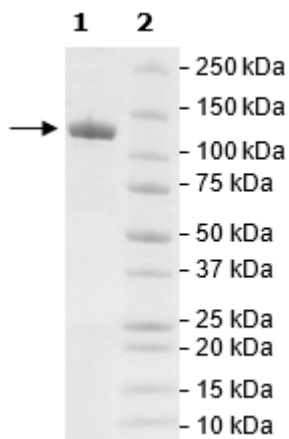


Product Information

Construct:	PARP1 (GST-2-1014(end))
Concentration:	1.20 mg/ml
Species:	Human
Formulated In:	40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.5 mM Glutathione, 3 mM DTT and 20% glycerol
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_001618
MW:	139 kDa
Purity:	≥90%
Assay Conditions:	Assay was performed according to PARP1 Chemiluminescent Assay Kit (BPS Bioscience #80551) with PARP1 serially diluted from 200 ng- 0 ng.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



PARP1 Activity

