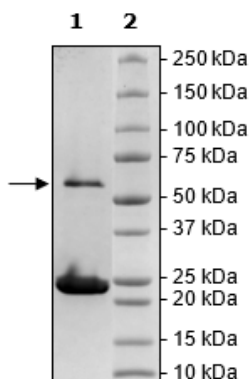


Product Information

Construct:	PARP11 (GST-8-338(end)-His)
Concentration:	0.21 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20 mM glutathione, 20% glycerol, and 3 mM DTT.
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_020367
MW:	65 kDa
Purity:	22%
Assay Conditions:	Assay done according to PARP11 Chemiluminescent Assay Kit (BPS Bioscience #80561) with various concentrations of PARP11.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PARP11 Activity

