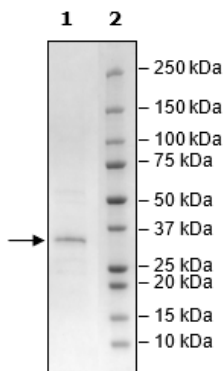


Product Information

Construct:	PARP7 (FLAG-400-657(end))
Concentration:	0.39 mg/ml
Species:	Human
Formulated In:	50 mM Tris, pH 8.0, 150 mM NaCl, 0.8 M Arginine, and 10% 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_015508
MW:	31 kDa
Purity:	79%
Assay Conditions:	Assay was performed according to PARP7 Chemiluminescent Assay Kit (BPS Bioscience #79729) with PARP7 titrated 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



PARP7 Activity

