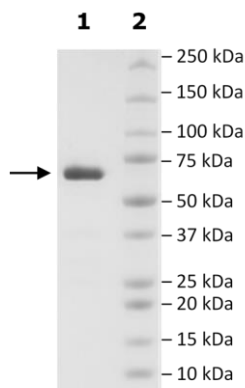


Product Information

Construct:	PARP14 (His-GST-TEV-1470-1801(end))
Concentration:	3.46 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, 200 mM imidazole
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_017554
MW:	67 kDa
Purity:	≥90%
Assay Conditions:	Assay was performed according to the PARP14 Chemiluminescent Assay Kit (BPS Bioscience #80568) with PARP14 serially titrated 2-fold from 1200 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



PARP14 Activity

