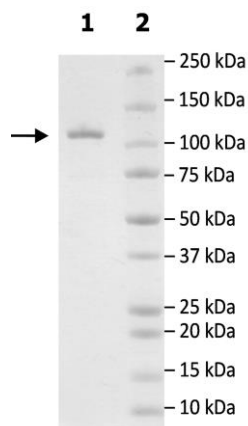


Product Information

Description:	Recombinant human PARP4 (poly(ADP-ribose) polymerase family member 4), encompassing amino acids 2-700. This construct contains an N-terminal His-tag (6xHis) followed by a GST-tag and a Thrombin Cleavage sight to facilitate removal of the GST-tag. This recombinant protein was affinity purified.
Species	Human
Construct:	PARP4 (His-GST-Th-2-700)
Concentration:	0.78 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT
MW:	107 kDa
Genbank Accession:	NM_006437
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was done according to PARP12 Chemiluminescent Assay Kit (BPS Bioscience #78504) with various amounts of PARP4.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PARP4 Activity

