

Product Information

Description:	Recombinant human PARP 4 (poly(ADP-ribose) polymerase family member 4), encompassing amino acids 2-700. This construct contains an N-terminal FLAG-Tag. This recombinant protein was affinity purified.
Species	Human
Construct:	PARP4 (FLAG-2-700)
Concentration:	0.67 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, 3mM DTT, and 100 µg FLAG peptide.
MW:	80 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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

