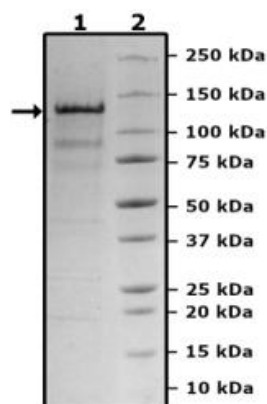


## Product Information

<b>Description:</b>	Recombinant human PARG (Poly(ADP-ribose) glycohydrolase), full length encompassing amino acids 2-976 (end). The construct contains an N-terminal His-Tag (6xHis). The recombinant protein was affinity purified.
<b>Background:</b>	PARG is the major enzyme responsible for the catabolism of poly(ADP-ribose), a reversible covalent-modifier of chromosomal proteins. It reverses PARP-mediated ADP-ribosylation and plays a role in DNA damage responses. PARG inhibition may prove beneficial in cancer therapy.
<b>Species:</b>	Human
<b>Construct:</b>	PARG (His-2-976(end))
<b>Concentration:</b>	0.40 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	70%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 610 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 0.5 mM TCEP, and 25 mM imidazole
<b>MW:</b>	112 kDa
<b>Genbank Accession:</b>	NM_003631.5
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Assay Conditions:</b>	Various concentrations of PARG were assayed using Fluorogenic Assay Kit (BPS Bioscience #78858) at room temperature with 4 μM TFMU-ADPr (substrate). The fluorogenic product was measured after 1 hour (Excitation: 385 nm / Emission: 502 nm). Enzyme concentration for IC <sub>50</sub> : 2.5 ng/rxn (~0.9 nM).
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PARG Activity

