

## PARG (aa 448-976)

**CATALOG NO.:** MSC-11-584

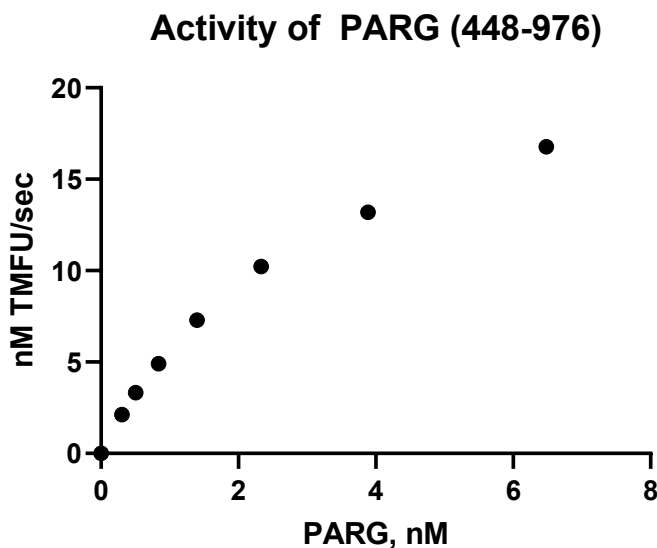
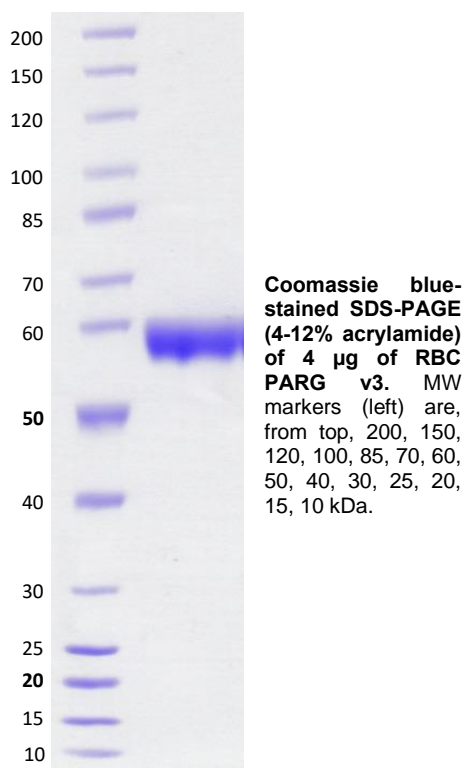
**LOT NO.:**

**DESCRIPTION:** Human recombinant PARG (residues 448-976; Genbank Accession # NM\_001303486; MW = 62.7 kDa) expressed as a C-terminal 8xHis-tagged protein in *E. coli*.

**PURITY:** >95% by SDS-PAGE

**SUPPLIED AS:** \_\_\_ µg/µL in 50 mM Tris, pH 8.0, 500 mM NaCl, 10% glycerol, 0.25 mM TCEP as determined by OD595.

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Activity of PARG (448-976) was measured in the presence of 50 µM TMFU-ADPr. Product formation was monitored at exc/em 385/502 nm using ClarioStar (BMG Labtech).

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com