

# TREVIGEN<sup>®</sup> Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

---

## Poly(ADP-ribose) glycohydrolase (PARG)

**Catalog #:** 4680-096-01

**Size:** 300  $\mu$ l

**Concentration:** 1  $\mu$ g/ml

**Description:** Poly(ADP-ribose) glycohydrolase (PARG) degrades poly(ADP-ribose) (PAR) polymers synthesized by poly(ADP-ribose) polymerase (PARP-1). PARG is ideal for use as a positive control in Trevigen's PARG Assay Kits (Cat# 4682-096-K and 4683-096-K) and in Western blot analysis of PARG in cell extracts.

**Source:** Purified from *E. coli* containing a recombinant plasmid harboring the catalytic domain of the bovine PARG gene.

**Storage Buffer:** 20 mM KPO<sub>4</sub> (pH 7.2), 50 mM KCl, 0.1 mg/ml BSA, 0.1% Triton<sup>®</sup> X-100, 1 mM DTT, 50% Glycerol.

**Measurement of Activity:** PARG activity is measured by the loss of biotinylated PAR from histones attached to the wells of a 96 well plate.

**Storage Conditions:** Stable for at least one year when stored at -20 °C.

### References:

1. Koh DW, Dawson VL, Dawson TM. 2005. The road to survival goes through PARG. *Cell Cycle*. 4:397-399.
2. Oei SL, Keil C, Ziegler M. 2005. Poly(ADP-ribosylation) and genomic stability. *Biochem Cell Biol*. 83:263-269.
3. Cuzzocrea S, Wang ZQ. 2005. Role of poly(ADP-ribose) glycohydrolase (PARG) in shock, ischemia and reperfusion. *Pharmacol Res*. 52:100-108.
4. Bonicalzi ME, Haince JF, Droit A, Poirier GG. 2005. Regulation of poly(ADP-ribose) metabolism by poly(ADP-ribose) glycohydrolase: where and when? *Cell Mol Life Sci*. 62:739-750.
5. Patel NS, Cortes U, Di Poala R, Mazzon E, Mota-Filipe H, Cuzzocrea S, Wang ZQ, Thiemermann C. 2005. Mice lacking the 110-kD isoform of poly(ADP-ribose) glycohydrolase are protected against renal ischemia/reperfusion injury. *J Am Soc Nephrol*. 16:712-719.
6. Patel CN, Koh DW, Jacobson MK, Oliveira MA. 2005. Identification of three critical acidic residues of poly(ADP-ribose) glycohydrolase involved in catalysis: determining the PARG catalytic domain. *Biochem J*. 388:493-500.

## TREVIGEN<sup>®</sup>

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: [info@trevigen.com](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

## Related Products Available From Trevigen

| Catalog #   | Description  | Size        |
|-------------|--|-------------|
| 4682-096-K  | HT Chemiluminescent PARG Assay Kit   | 96 tests    |
| 4683-096-K  | HT Colorimetric PARG Assay Kit   | 96 tests    |
| 4520-096-K  | PARP in vivo Pharmacodynamic Assay II  | 96 tests    |
| 4676-096-K  | HT Universal Chemiluminescent PARP Assay Kit/w<br>Histone Coated Strip Wells | 96 Samples  |
| 4677-096-K  | HT Universal Colorimetric PARP Assay Kit/w Histone Coated<br>Strip Wells     | 96 Samples  |
| 4684-096-K  | HT Colorimetric PARP Apoptosis Assay   | 96 tests    |
| 4685-096-K  | HT Chemiluminescent PARP Apoptosis Assay                                     | 96 tests    |
| 4690-096-K  | HT F Homogeneous PARP Inhibition Assay Kit                                   | 96 tests    |
| 4668-100-01 | Recombinant Human PARP (High Specific Activity)                              | 1000 Units  |
| 4667-250-EB | Recombinant Human PARP Enzyme  | 250 $\mu$ l |
| 4667-50-03  | 3-Aminobenzamide PARP inhibitor (200 mM)                                     | 60 $\mu$ l  |
| 4667-50-11  | Benzamide PARP inhibitor (8 mM)  | 100 $\mu$ l |
| 4667-50-10  | 6(SH)-Phenanthridinone PARP inhibitor (160 $\mu$ M)                          | 100 $\mu$ l |
| 4667-50-9   | 4-Amino-1,8-naphthalimide PARP inhibitor (800 $\mu$ M)                       | 100 $\mu$ l |

### PARG

Catalog#: 4680-096-01

Storage: -20 °C

**TREVIGEN®**  
1-800-873-8443