

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Poly(ADP-ribose) glycohydrolase (PARG)

Catalog #: 4680-096-01

Size: 300 µl
Concentration: 1 µg/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 KPO₄ (pH 7.2), 50 mM KCl, 0.1 mg/ml BSA, 0.1% Triton® 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.

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 PARG Assay Kit/w Histone Coated Strip Wells	96 Samples
4677-096-K	HT Universal Colorimetric PARG Assay Kit/w Histone Coated Strip Wells	96 Samples
4684-096-K	HT Colorimetric PARG Apoptosis Assay	96 tests
4685-096-K	HT Chemiluminescent PARG Apoptosis Assay	96 tests
4690-096-K	HT F Homogeneous PARG Inhibition Assay Kit	96 tests
4668-100-01	Recombinant Human PARG (High Specific Activity)	1000 Units
4667-250-EB	Recombinant Human PARG Enzyme	250 µl
4667-50-03	3-Aminobenzamide PARG inhibitor (200 mM)	60 µl
4667-50-11	Benzamide PARG inhibitor (8 mM)	100 µl
4667-50-10	6(SH)-Phenanthridinone PARG inhibitor (160 µM)	100 µl
4667-50-9	4-Amino-1,8-naphthalimide PARG inhibitor (800 µM)	100 µl

© 2010 Trevigen, Inc. All Rights Reserved. Triton is a registered trademark of Union Carbide Chemicals and Plastic Company. Trevigen is a registered trademark of Trevigen, Inc. E8/26/10v1

TREVIGEN®

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 • www.trevigen.com

PARG

Catalog#: 4680-096-01

Storage: -20 °C

TREVIGEN®
1-800-873-8443