

# TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## Human Poly(ADP-Ribose) Polymerase (PARP)

**Catalog #:** 4667-250-EB

**Contents:** 4667-250-01 PARP-1 **Size:** 250 µl  
4671-096-02 20X PARP Buffer 500 µl

**Description:** Poly ADP-ribosylation by PARP of nuclear proteins is a posttranslational event that occurs in response to DNA damage. The enzyme provided is less than 50% pure by SDS-PAGE and Coomassie staining. The enzyme is useful as a positive control for both Western blot analysis of ribosylated proteins and PARP activity assays.

**Source:** Purified from *E. coli* containing a recombinant plasmid harboring the human PARP-1 gene.

**Unit Definition:** One unit of PARP-1 incorporates 10 femtomoles of NAD onto immobilized histones in 30 minutes. Number of units provided (>5 Units/ µl) is lot specific.

**Storage Buffer:** 20 mM Tris-Cl, pH 8.0, 200 mM NaCl, 1 mM DTT, and 50% glycerol.

**Storage Conditions:** Store at -20°C in a manual defrost freezer. For long term storage, freeze in working aliquots at -80°C. Avoid freeze thaw cycles.

**References:** 1. Satoh, M.S. and T. Lindahl. 1994. Role of poly(ADP-ribose) formation in DNA repair. *Nature* 356:356-358.  
2. Lautier, D., J. Lagueux, J. Thibodeau, L. Menard, and G.G. Poirier. 1993. Molecular and biochemical features of poly(ADP-ribose) metabolism. *Mol Cell Biochem* 122:171-193.

**Lot Specific Data:** see reverse

## Related Products Available From Trevigen

Catalog #	Description	Size
4668-100-01	Recombinant Human PARP (High Specific Activity)	1000 Units
4667-50-EB	Recombinant Human PARP Enzyme	50 µl
4520-096-K	PARP in vivo Pharmacodynamic Assay II	96 tests
4676-096-K	HT Universal Chemiluminescent PARP Assay Kit/w Histone Coated Strip Wells	96 Samples
4677-096-K	HT Universal Colorimetric PARP Assay Kit/w Histone Coated 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
4682-096-K	HT Chemiluminescent PARG Assay Kit	96 tests
4683-096-K	HT Colorimetric PARG Assay Kit	96 tests
4667-50-03	3-Aminobenzamide PARP inhibitor (200 mM)	60 µl
4667-50-11	Benzamide PARP inhibitor (8 mM)	100 µl
4667-50-10	6(SH)-Phenanthridinone PARP inhibitor (160 µM)	100 µl
4667-50-9	4-Amino-1,8-naphthalimide PARP inhibitor (800 µM)	100 µl

**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

**Human Poly(ADP-ribose) Polymerase (PARP-1)**

**Catalog #: 4667-250-EB**

**Storage: -80 °C**

**TREVIGEN®**  
**1-800-873-8443**