

# TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

## Human Poly(ADP-Ribose) Polymerase (PARP): High Specific Activity

Catalog #: 4668-100-01

Size: 1,000 Units

**Description:** Poly ADP-ribosylation by PARP of nuclear proteins is a posttranslational event that occurs in response to DNA damage. PARP inhibition prevents tissue damage in animal models of myocardial and neuronal ischemia, diabetes, septic shock, and vascular stroke. High specific activity PARP is ideal for use as a positive control in Trevigen's PARP Assay Kits (e.g. cat#s 4677-096-K; 4684-096-K; 4690-096-K) and in Western blot analysis of PARP in cell extracts.

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

**Unit Definition:** One unit of PARP incorporates 10 femtomoles of NAD onto 5 µg of immobilized histones in 30 minutes.

**Assay Conditions:** 50 mM Tris-Cl (pH 8.0), 25 mM MgCl<sub>2</sub>, 34 µM of a mixture of NAD and biotinylated NAD, 0.02 mg/ml activated DNA, and serial dilutions of PARP in a reaction volume of 50 µl. Incubation is for 30 minutes at 22 °C. The incorporation of biotinylated poly(ADP-ribose) is quantified by incubation with Streptavidin-HRP followed by addition of a colorimetric HRP substrate and measurement of the absorbance at 450 nm.

**Storage Buffer:** 20 mM Tris-Cl (pH 8.0), 200 mM NaCl, 1 mM DTT, 0.1% Triton X-100, 50% glycerol, and 0.1 mg/ml BSA.

**Storage Conditions:** The enzyme retains over 66% of its activity after 24 hr at 37 °C. Stable for at least one year when stored at -20 °C.

### 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.
3. Bazar, L.S. Quantitation of PARP Activity in cell extracts and determination of PARP inhibitor efficacy. *Bioscience Technology*, July 2005, 44-48.
4. Curtin NJ. 2005. PARP inhibitors for cancer therapy. *Expert Rev Mol Med*. **7**:1-20.
5. Kim MY, Mauro S, Gevry N, Lis JT, Kraus WL. 2004. NAD-Dependent Modulation of Chromatin Structure and Transcription by Nucleosome Binding Properties of PARP-1. *Cell* **119**:803-814.
6. Virag, L., and Szabo, C. 2002. The therapeutic potential of Poly(ADP-Ribose) Polymerase inhibitors. *Pharmacological Reviews* **54**:375-429.

**Lot Specific Data:** see reverse

© 2008 Trevigen, Inc. All Rights Reserved. Trevigen is a registered trademark of Trevigen, Inc.

E8/25/08v1

**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

Lot Specific Data:

Lot #:

Activity:

Protein Concentration:

Specific Activity:

**Human PARP  
High Specific Activity**

Catalog #: 4668-100-01

Storage: -20 °C

**TREVIGEN®  
1-800-873-8443**