

# TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

---

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

<b>Catalog #:</b>	4667-50-01	PARP-1	<b>Size:</b> 50 µl
	4667-250-01	PARP-1	250 µl
	4667-50-02	10X PARP Buffer	700 µl

**Description:** The enzyme is useful as a positive control for both Western blot analysis of ribosylated proteins and the study of PARP inhibitors.

**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 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, 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

## 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](mailto:info@trevigen.com) • [www.trevigen.com](http://www.trevigen.com)

Size:

Lot Specific Data:

Activity:

[Protein]:

Specific Activity:

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

**Catalog #: 4667-250-01**

**Lot #: 11611J6**

**Storage: -80 °C**

**TREVIGEN®**  
**1-800-873-8443**