

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures.

Human Poly(ADP-Ribose) Polymerase (PARP)

Catalog #: 4667-50-01 PARP

4667-50-02 10X PARP Buffer

Size: 50 µl

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 gene.

Unit Definition: One unit of PARP incorporates 100 pmoles of poly(ADP) from NAD into acid-insoluble form in 1 minute at 22°C.

Activity: Greater than or equal to 2.0 Units/µl.

Assay Conditions: 1X PARP Buffer (50 mM Tris-Cl (pH 8.0), and 25 mM MgCl₂), 1 mM ³²P-NAD, 0.1 mg/ml activated salmon testes DNA, 0.1 mg/ml Histone H1, and 1 µl of enzyme in a reaction volume of 100 µl. After incubating for 1 minute, 5 minutes, and 10 minutes at 22°C, the reactions are stopped by the addition of 1 ml of 20% trichloroacetic acid. Acid-insoluble counts are then measured for quantitation of unit activity.

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

Storage Conditions: Store at -20°C. For long term storage, freeze in working aliquots at -80°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.

Lot Specific Data: see reverse

Lot Specific Data:

Lot #: enquire for lot specific information

Activity:

Total Protein:

Specific Activity:

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

Catalog #: 4667-50-01

Storage: -20°C

TREVIGEN[®]

1-800-873-8443

TREVIGEN[®]

8405 Helgerman Court, Gaithersburg, MD 20877 USA

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

Fax: 301-216-2801 • e-mail: info@trevigen.com • www.trevigen.com