

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Poly(ADP-ribose) (PAR) Polymer

Catalog #: 4336-100-01

Volume: 100 µl

Concentration: 10 µM

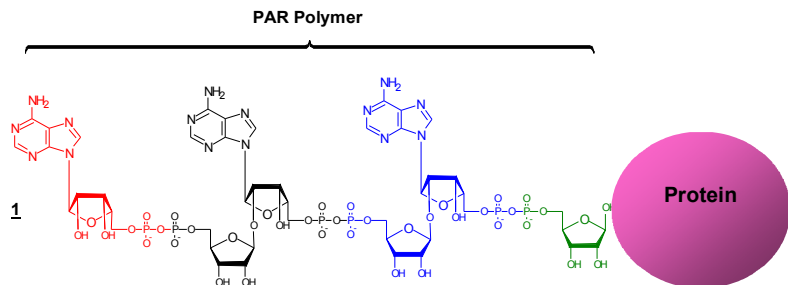


Figure 1: Poly (ADP-ribose) (PAR)

Description: PAR (figure 1) was synthesized using poly(ADP-ribose) polymerase (PARP) in the presence of NAD, and purified by gel filtration. The PAR chain length ranges from 2 to 300 monomers, and it is recognized by Trevigen's anti-PAR polymer monoclonal antibody (Cat# 4335-MC-100) in a PAR ELISA (figure 2). PAR concentration was determined using the following equation:¹

$$[\text{PAR}] = \frac{(A_{258}) \text{ cm}^{-1}}{13,500 \text{ cm}^{-1} \text{ M}^{-1}}$$

Storage Conditions: PAR is provided in 10 mM Tris-HCl (pH 8.0), 1 mM EDTA and is stable for at least 1 year at -80 °C. It may be aliquoted to avoid repeated freeze-thawing.

Applications: Immunodetection as a standard in PAR ELISAs. Empirical determination will be required for optimal results.

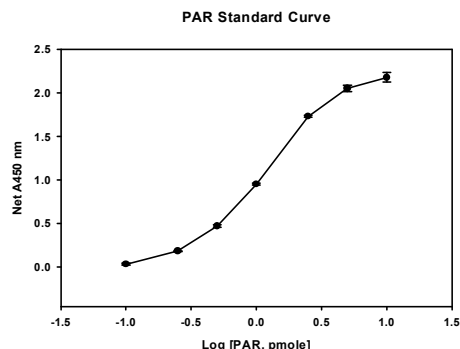


Figure 2: Serial dilutions of PAR were applied in triplicate to an anti-PAR pre-coated 96-well plate and following incubation for 1.5 hrs, rabbit polyclonal anti-PAR (Cat# 4336-BPC-100) was added and the plate was incubated at room temperature for 1.5 hrs. After washing, goat anti-rabbit IgG-HRP was added and incubated at room temperature for 1 hr. Post washing, wells were developed using TACS-sapphire™ (Cat# 4822-96-08) for 5 minutes, and the reaction was stopped using phosphoric acid. The A₄₅₀ was determined using a plate reader.

© 2009 Trevigen, Inc. All Rights Reserved. Trevigen is a registered trademark of Trevigen, Inc.

E9/11/09v1

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

References:

1. Shah, G.M., et al. 1995. Methods for biochemical study of poly(ADP-ribose) metabolism *in vitro* and *in vivo*. *Anal Biochem* **227**:1-13.
2. Affar, E.B., et al. 1998. Immunodot blot method for the detection of poly(ADP-ribose) synthesized *in vitro* and *in vivo*. *Anal Biochem* **259**:280-283.
3. Menard, L. and G.G. Poirier. 1987. Rapid assay of poly(ADP-ribose) glycohydrolase. *Biochem Cell Biol* **65**:668-673.

**Poly(ADP-ribose)
(PAR) polymer**

Catalog#: 4336-100-01

Storage: -80 °C

TREVIGEN®
1-800-873-8443