

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Biotin (Terminal) – Poly ADP-ribose (PAR) Polymer

Catalog #: 4336-100-02

Volume: 100 µl

Concentration: 10 µM

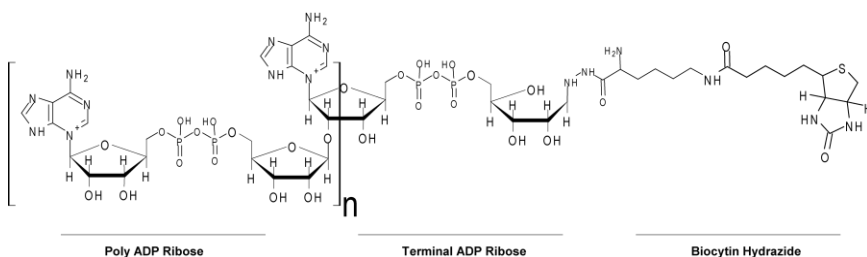


Figure 1: Terminally biotinylated Poly ADP-ribose (PAR)

Description: Biotin (Terminal)-PAR (figure 1) was synthesized by reductive amination¹ of pure PAR polymer. Free biotin was removed by desalting over Sephadex G25 column. The PAR chain length³ ranges from 2-17 mer² (28%), 17-34 mer (34%), > 34 mer (38%), and it is recognized by Trevigen's PAR polymer monoclonal antibody (cat# 4335-MC-100) and polyclonal antibodies (cat# 4336-BPC-100). The Biotin (Terminal)-PAR concentration was determined using the following equation⁴.

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

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

Applications: Co-precipitation assay, ELISA and Western blot analysis.

References:

1. Fahrner, J., Kranaster, R., Altmeyer, M., Marx, A., Bürkle, A., 2007. Quantitative analysis of the binding affinity of poly(ADP-ribose) to specific binding proteins as a function of chain length. *Nucleic Acids Res* 35, e143.
2. As the material was desalted over a Sephadex G25 column, small oligomers of ADP ribose (2-10 mer), will be absent in the final product.
3. Tan, Edwin S, Kristin A Krukenberg, and Timothy J Mitchison. 2012. "Large-scale Preparation and Characterization of poly(ADP-ribose) and Defined Length Polymers." *Analytical Biochemistry* **428**: 126–136.
4. 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.

© 2012 Trevigen, Inc. All Rights Reserved. Trevigen is a registered trademark of Trevigen, Inc.

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

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

Catalog#: 4336-100-01

Storage: -80 °C

TREVIGEN®

1-800-873-8443