

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Anti-PAR Monoclonal Affinity Purified

Catalog #: 4335-AMC-050

Size: 50 µl

Description: Trevigen's affinity purified mouse monoclonal antibody (cat# 4335-AMC-050) is specific for poly (ADP-ribose) (PAR) polymer. It can be used to detect ribosylated proteins by ELISA, Western blot, immunocytochemistry *in situ*, and for immunopurification.

Storage Conditions: Store the anti-PAR monoclonal at -20°C (manual defrost freezer).

Physical State: This antibody is an affinity purified immunoglobulin from mouse ascites in 1X PBS containing 50% glycerol at 1 mg/ml.

Immunogen: Purified ADP-ribose polymers between 2 and 50 units long.

Ig Class: IgG₃

Specificity: The antibody is specific for PAR polymers 2 to 50 units long, but does not recognize structurally related RNA, DNA, ADP-ribose monomers, NAD, or other nucleic acid monomers.

Applications: ELISA, Western analysis, immunoprecipitation, and immunopurification. For Western blots, an antibody dilution of 1:1000 is recommended.

References:

1. Kupper, J.H., G. de Murcia, and A. Burkle. 1990. Inhibition of poly(ADP-ribosylation) by overexpressing the poly(ADP-ribose) polymerase DNA binding domain in mammalian cells. *J Biol Chem* **265**: 18721-18724.
2. Malik, N., M. Miwa, T. Sugimura, P. Thraves, and M. Smulson. 1983. Immunoaffinity fractionation of the poly(ADP-ribosyl)ated domains of chromatin. *PNAS USA* **80**:2554-2558.
3. Smulson, M.E. et al.. 1983. In ADP-Ribosylation, DNA Repair and Cancer (Miwa, M., O. Hayaishi, S. Shall, M. Smulson, and T. Sugimura, eds.), pp. 49-70. Tokyo:Japan *Scientific Societies Press*.
4. Kawamitsu, H., H. Hoshino, H. Okada, M. Miwa, H. Momoi, and T. Sugimura. 1984. Monoclonal antibodies to poly(adenosine diphosphate ribose) recognize different structures. *Biochem* **23**:3771- 3777.
5. 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.

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

**Anti-PAR
Monoclonal
Affinity Purified
Cat#: 4335-AMC-050
Storage: -20°C
(Manual Defrost Freezer)**
TREVIGEN®
1-800-873-8443