



TREVIGEN®

Products for PARP/PARG Research

Poly ADP-ribosylation of nuclear proteins is a post-translational event that occurs in response to DNA damage. Poly-ADP-ribose polymerase (PARP) catalyzes the NAD-dependent addition of poly-ADP-ribose (PAR) to adjacent nuclear proteins. PARP plays an important role in DNA repair but can also lead to apoptosis by depleting the cellular NAD pool. PARP inhibition has been shown to prevent tissue damage in animal models of myocardial and neuronal ischemia, diabetes, septic shock, and vascular stroke.

Standard and High Specific Activity PARP Enzyme

Trevigen offers a “standard” PARP purified to ~50% purity and a high specific activity (HSA) enzyme purified to greater than 95%. The standard PARP enzyme is useful as a positive control for Western blot analysis of ribosylated proteins and PARP Activity Assays. HSA PARP is recommended for detailed studies and drug discovery applications.

Description	Size	Catalog #
Recombinant Human PARP Enzyme	250 µl	4667-250-01
Recombinant Human PARP Enzyme	1 ml	4667-1K-01
High-Specific Activity Recombinant Human PARP Enzyme	1,000 units	4668-100-01
High-Specific Activity Recombinant Human PARP Enzyme	5,000 units	4668-500-01
High-Specific Activity Recombinant Human PARP Enzyme	20,000 units	4668-2K-01

PARP and PAR Antibodies

The PARP monoclonal antibody can be used for analysis of PARP protein levels and cleavage of PARP in apoptotic cells. It recognizes both the full length protein (116 kDa) and an 85 kDa cleavage fragment associated with apoptosis. The PAR antibody is specific for PAR polymers synthesized by a variety of PARP isoforms (PARP-1,-2,-3, tankyrase, vPARP, and others).

Description	Size	Catalog #
PARP MAb (Clone C2-10)	50 µg	4338-MC-50
PAR MAb (Clone 10HA)	100 µl	4335-MC-100
PAR MAb (Clone 10HA)	1000 µl	4335-MC-1K
Poly(ADP) Ribose (PAR) Polymer	100 µl	4336-100-01
PAR PAb (Rabbit)	100 µl	4336-BPC-100

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS



TREVIGEN®

ANGIOGENESIS • DNA DAMAGE & REPAIR • APOPTOSIS

Products for PARP/PARG Research

PARP and PARG Kits

The PARP Assay Kits measure incorporation of biotinylated Poly (ADP-ribose) onto histone proteins in 96-well plates. The assays allow screening of PARP inhibitors and measurement of PARP activity in cell and tissue extracts. PARG Assay Kits allow screening of PARG inhibitors and measurement of PARG activity in cell extracts.

Description	Size	Detection	Sensitivity	Catalog #
PARP Assay Kit	1 plate	Colorimetric	1 m unit	4677-096-K
PARP Assay Kit	96 Tests	Colorimetric	1 m unit	4671-096-K
PARP Assay Kit	1 plate	Chemiluminescent	2.5 m units	4676-096-K
PARP Assay Kit	96 Tests	Chemiluminescent	2.5 m units	4675-096-K
HT PARP Apoptosis Assay Kit	96 Tests	Colorimetric	0.1 m unit	4684-096-K
HT PARP Apoptosis Assay Kit	1 plate	Chemiluminescent	0.1 m unit	4685-096-K
PARG Assay Kit	1 plate	Colorimetric	0.5 pg	4683-096-K
PARG Assay Kit	1 plate	Chemiluminescent	0.5 pg	4682-096-K
Histone-coated Strip Well Plate (natural)	1 plate	N/A	N/A	4677-096-P
Histone-coated Strip Well Plate (white)	1 plate	N/A	N/A	4678-096-P