

PARP-1 N-terminal antibody (pAb)

Catalog Nos: 39559, 39061, 39560

RRID: AB_2793257

Isotype: Serum

Application(s): ChIP, WB

Reactivity: Human, Mouse

Volumes: 100 μ l, 50 μ l, 10 μ l

Purification: None

Host: Rabbit

Molecular Weight: 120 kDa

Background: PARP-1 N-terminal (ADPRT) encodes a chromatin-associated enzyme, poly(ADP-ribose)transferase, that modifies various nuclear proteins by poly(ADP-ribose)ation. The modification is dependent on DNA and is involved in the regulation of various important cellular processes such as differentiation, proliferation and tumor transformation. It also plays a role in the regulation of the molecular events involved in the recovery of cells from DNA damage. Cleavage of PARP-1 (ADPRT) occurs following caspase activation during apoptosis.

For additional information on PARP-1, please see the review article PARP-1: An Abundant and Ubiquitous Protein with Roles in Many Cellular Processes in the Targets & Applications section or our website.

Immunogen: This PARP-1 N-terminal antibody was raised against a His-Tagged fusion protein corresponding to the N-terminal half of human PARP-1 was used to generate this PARP-1 antibody.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic. For your convenience, an IgG version (Catalog No. 61639) of this antibody that was purified by Protein A Chromatography is also available.

Application Notes:

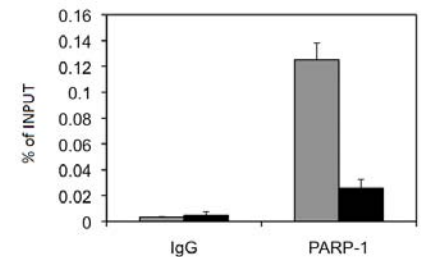
Applications Validated by Active Motif:

ChIP: 8 μ l per ChIP

WB: 1:5,000 - 1:50,000 dilution

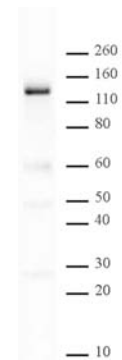
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



PARP-1 N-terminal antibody tested by ChIP analysis.

Chromatin IP performed with chromatin from 3T3 cells using 8 μ l of PARP-1 N-terminal antibody or rabbit IgG as a negative control. Real time, quantitative PCR (RT-qPCR) was performed on DNA purified from each of the ChIP reactions using a primer pair specific for the PPARG2 gene. Data are presented as percent of input from cells with a PARP-1 knock-down (black bars) or a control knock-



PARP-1 N-terminal antibody tested by Western blot.

HeLa cell nuclear extract (20 μ g per lane) was probed with PARP-1 N-terminal antibody at a dilution of 1:20,000.