

PARP-1 C-terminal antibody (pAb)

Catalog Nos: 39561, 39562

RRID: AB_2793258

Application(s): ChIP, ChIP-Seq, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: None

Host: Rabbit

Isotype: Serum

Molecular Weight: 120 kDa

Background: PARP-1 C-terminal (ADPRT) encodes a chromatin-associated enzyme, poly(ADP-ribosyl)transferase, that modifies various nuclear proteins by poly(ADP-ribosyl)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 C-terminal antibody was raised against a His-Tagged fusion protein corresponding to the C-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.

Application Notes:

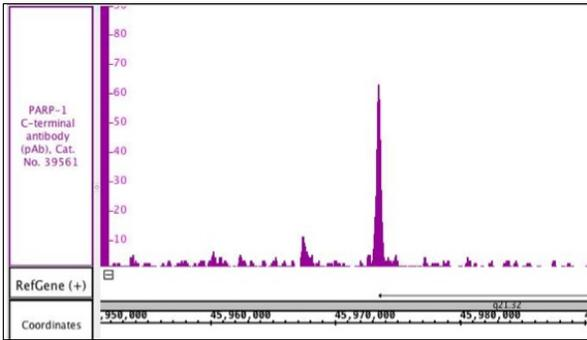
Applications Validated by Active Motif:

ChIP: 10 µl per ChIP

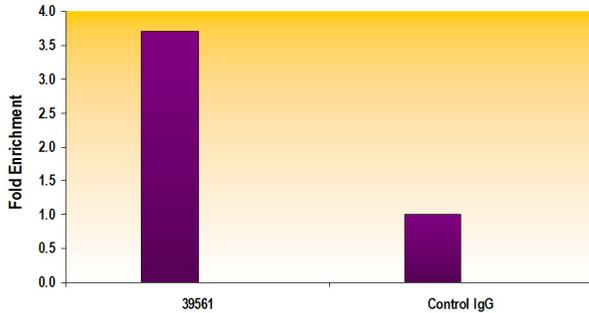
WB: 1:1,000 - 1:6,000 dilution

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.

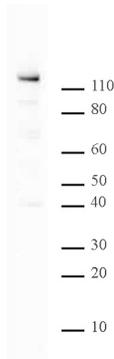


PARP-1 C-terminal antibody (pAb) antibody (rAb) tested by ChIP-Seq. Chromatin immunoprecipitation (ChIP) was performed using the **ChIP-IT[®] High Sensitivity Kit** (Cat. No. 53040) with 30 µg of chromatin from human NCI-H209 lung cancer cells and 10 µl of PARP-1 C-terminal antibody (pAb). ChIP DNA was sequenced on the Illumina NextSeq and 7.7 million sequence tags were mapped to identify PARP binding sites on chromosome 17.



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

Chromatin IP performed using the ChIP-IT[®] Express Kit (Catalog No. 53008) and MCF-7 chromatin (1.5 x 10⁶ cell equivalents per ChIP) using 10 µl of PARP-1 C-terminal antibody or the equivalent amount of 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 BRCA2 gene promoter. Data are presented as Fold Enrichment of the ChIP antibody signal versus the negative control IgG using the ddCT method.



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

HeLa nuclear extract (20 µg per lane) was probed with PARP-1 C-terminal antibody at a dilution of 1:6,000.