

## CDKN2A antibody (pAb)

**Catalog Nos:** 61619, 61620

**RRID:** AB\_2793705

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100  $\mu$ l, 10  $\mu$ l

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 16 kDa

**Background:** CDKN2A (Cyclin-Dependent Kinase Inhibitor 2A) is capable of inducing cell cycle arrest in G1 and G2 phases. Acts as a tumor suppressor. Binds to MDM2 and blocks its nucleocytoplasmic shuttling by sequestering it in the nucleolus. This inhibits the oncogenic action of MDM2 by blocking MDM2-induced degradation of p53 and enhancing p53-dependent transactivation and apoptosis. Also induces G2 arrest and apoptosis in a p53-independent manner by preventing the activation of cyclin B1/CDC2 complexes. Binds to BCL6 and down-regulates BCL6-induced transcriptional repression. Binds to E2F1 and MYC and blocks their transcriptional activator activity but has no effect on MYC transcriptional repression. Binds to TOP1/TOPOI and stimulates its activity. This complex binds to rRNA gene promoters and may play a role in rRNA transcription and/or maturation. Interacts with NPM1/B23 and promotes its polyubiquitination and degradation, thus inhibiting rRNA processing. Interacts with COMMD1 and promotes its 'Lys63'-linked polyubiquitination. Interacts with UBE2I/UBC9 and enhances sumoylation of a number of its binding partners including MDM2 and E2F1. Binds to HUWE1 and represses its ubiquitin ligase activity. May play a role in controlling cell proliferation and apoptosis during mammary gland development. Isoform 6 may be involved in regulation of autophagy and caspase-independent cell death; the short-lived mitochondrial isoform is stabilized by C1QBP.

Multiple protein isoforms are encoded by this gene. The p16 isoform, also known as INK4a, is a CDK4 & CDK6 kinase inhibitor. It functions as negative regulator of the proliferation of normal cells and plays a role in aging and senescence.

**Immunogen:** This antibody was raised against a peptide within the N-terminal region of the p16/INK4a isoform of human CDKN2A.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

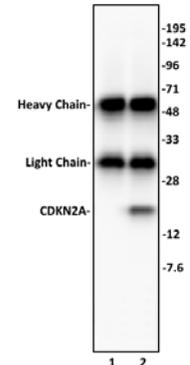
### Application Notes:

Applications Validated by Active Motif:

IP: 10  $\mu$ l per IP

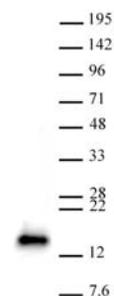
WB: 1:500 - 1:2,000 dilution

**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.



### CDKN2A antibody (pAb) tested by Immunoprecipitation.

10  $\mu$ l of CDKN2A antibody was used to immunoprecipitate CDKN2A from 400  $\mu$ g of HeLa whole cell extract (lane 2). 10  $\mu$ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the CDKN2A antibody at a dilution of 1:500.



### CDKN2A antibody (pAb) tested by Western blot.

Detection of CDKN2A by Western blot. The analysis was performed using 30  $\mu$ g HeLa whole-cell extract and CDKN2A (pAb) at a 1:500 dilution.