

## PP2A antibody (pAb)

**Catalog Nos:** 39192, 39193

**RRID:** AB\_2793178

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

**Reactivity:** Human, Mouse, Rat, Wide Range Predicted

**Volumes:** 200 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Isotype:** IgG

**Molecular Weight:** 36 kDa

**Background:** PP2A – Protein Phosphatase 2A beta subunit is the catalytic subunit of protein phosphatase 2A. PP2A is involved in the regulation of a number of signaling pathways, and it is implicated in the negative control of cell growth and division. Mutations of PP2A occurs in many types of cancers.

**Immunogen:** This PP2A antibody was raised against a peptide corresponding to amino acids 288-303 of human PP2A.

**Buffer:** PBS containing 30% glycerol, 0.035% sodium azide and 1 mg/ml BSA. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

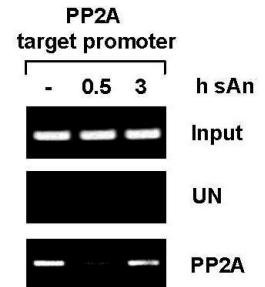
ChIP: 2 - 10 µl per ChIP

WB\*: 1:250 - 1:2,000 dilution>

\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

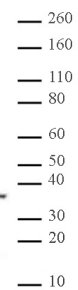
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This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



### PP2A antibody (pAb) tested by ChIP analysis.

Chromatin IP performed using Swiss 3T3 Chromatin (approximately 5 million cells treated with anisomycin and fixed at 0.5 and 3 hours) using 2.5 µl of PP2A antibody or the equivalent amount of IgG as a negative control.



### PP2A pAb tested by Western blot.

Nuclear extract of HeLa cells was probed with PP2A pAb (1:1,000 dilution).