

YAP1 antibody (pAb)

Catalog Nos: 61293, 61294

RRID: AB_2614990

Isotype: IgG

Application(s): IP, WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Concentration: 1.4 μ g/ μ l

Molecular Weight: 75 kDa

Background: YAP1 (Yes-associated protein 1, YAP, YKI, YAP2, YAP65) is a transcriptional regulator which can act both as a coactivator and a corepressor and is the critical downstream regulatory target in the Hippo signaling pathway that plays a pivotal role in organ size control and tumor suppression by restricting proliferation and promoting apoptosis. Is a critical modulator of epidermal stem cell proliferation. The core of this pathway is composed of a kinase cascade wherein MST1/MST2, in complex with its regulatory protein SAV1, phosphorylates and activates LATS1/2 in complex with its regulatory protein MOB1, which in turn phosphorylates and inactivates YAP1 oncoprotein and WWTR1/TAZ. Plays a key role to control cell proliferation in response to cell contact. Phosphorylation of YAP1 by LATS1/2 inhibits its translocation into the nucleus to regulate cellular genes important for cell proliferation, cell death, and cell migration. The presence of TEAD transcription factors are required for it to stimulate gene expression, cell growth, anchorage-independent growth, and epithelial mesenchymal transition (EMT) induction.

Immunogen: This YAP1 antibody was raised against a peptide corresponding to the C-terminal region of human YAP1.

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

Application Notes:

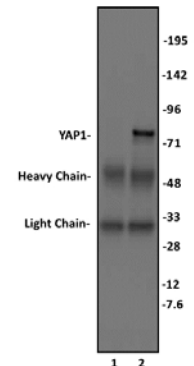
Applications Validated by Active Motif:

IP: 10 μ l per IP

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

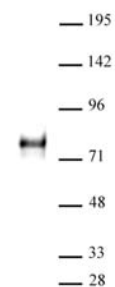
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



YAP1 antibody (pAb) tested by Immunoprecipitation.

10 μ l of YAP1 antibody was used to immunoprecipitate YAP1 from 400 μ g of HeLa whole cell extract (lane 2). 10 μ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the YAP1 antibody at a dilution of 1:500.



YAP1 antibody (pAb) tested by Western blot.

YAP1 detection by Western blot. The analysis was performed using 40 μ g HeLa whole-cell extract and YAP1 at a 1:500 dilution.