

p53 antibody (pAb)

Catalog Nos: 61657, 61658**RRID:** AB_2793723**Isotype:** IgG**Application(s):** WB**Reactivity:** Human**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 53 kDa

Background: p53 is the most important tumor suppressor in the genome. It is responsive to numerous genotoxic stresses, which activates its transcription factor activity, in turn causing cell-cycle arrest by activating expression of p21 Cip/WAF. Mutant p53 that has lost its DNA-binding function interferes with the activity of native p53 and leads to oncogenic transformation. Alternatively, transformation may be caused by overexpression of Mdm2/Hdm2, a ubiquitin ligase specific for p53, which causes its destabilization. Inactivation of p53 is often coincident with hyperactivation of NFκB (NFκB p50 and NFκB p65), both of which serve to inhibit apoptosis.

Immunogen: This antibody was raised against recombinant human p53 protein.

Buffer: Purified IgG in PBS with 50% glycerol and 0.02% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

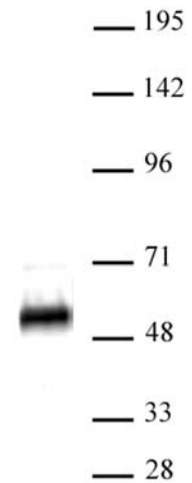
WB*: 1:500 - 1:2,000 dilution

*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.

For p53, we also offer AbFlex® p53 Recombinant Antibody (rAb). For details, see Catalog No. 91247.

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.

**p53 antibody (pAb) tested by Western blot.**

Detection of p53 by Western blot. HEK293 nuclear extract (30 µg) probed with p53 antibody at a 1:500 dilution.