

p53 antibody (mAb)

Catalog Nos: 39553, 39554

RRID: AB_2793254

Clone: DO1

Isotype: IgG

Application(s): ChIP, IP, WB

Reactivity: Human

Quantities: 200 µg, 10 µg

Purification: Protein G Chromatography

Host: Mouse

Concentration: 1 µg/µl

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 p53 antibody was raised against full-length recombinant human p53 protein.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

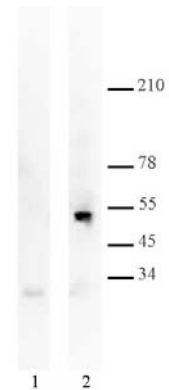
IP: 5 µg per IP

WB: 0.5 - 2 µg/ml dilution

This antibody is also available as an AbFlex® engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91247.

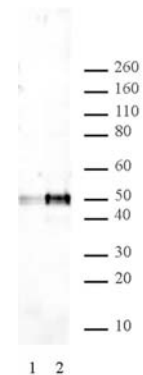
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product 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 mAb (Clone DO1) tested by Immunoprecipitation.

5 µg of p53 mAb (Clone DO1) was used to immunoprecipitate p53 from 500 µg of camptothecin-treated U2OS nuclear extract (lane 2). 5 µg of mouse IgG was also used as a control (lane 1). The immunoprecipitated protein was detected by western blotting using the same p53 antibody.



p53 mAb (Clone DO1) tested by Western blot.

Western Blot: Nuclear extract of U2OS cells (20 µg) probed with p53 mAb (Clone DO1) (1 µg/ml).

Lane 1: No treatment.

Lane 2: Cells treated with camptothecin.