

p53 antibody (mAb)

Mouse Anti-Human/Mouse/Rat p53

Instruction Manual

Catalog Number	PK-AB577-3036M
Description	p53, a 53 kDa protein, binds to a DNA consensus sequence, the p53 response element, and regulates normal cell cycle events by activating transcription of genes involved either in progression through the cell cycle, or causing arrest in G1 when the genome is damaged. In most transformed and tumor cells the concentration of p53 is increased 5-1000 fold over the concentration in normal cells, principally due to the increased half-life (4 hrs) compared to that of wild-type (20 min). p53 localizes in the nucleus, but is detectable at the plasma membrane during mitosis and certain mutations also modulate cytoplasmic/nuclear distribution. p53 downregulates Bcl-2 expression and upregulates Bax expression, but may not always be necessary for apoptosis.
Quantity	100 µg
Source / Host	Mouse
Immunogen	NA
Clone / IgG Subtype	Mouse IgG1; clone C6A5
Species Reactivity	human, mouse, rat
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) in PBS containing 1 mg/ml BSA and 1.5 mM sodium azide and 50% glycerol.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blotting (0.5-4 µg/ml), immunoprecipitation (4-8 µg/ml), Immunohistochemistry (4-8 µg/ml), immunocytochemistry (4-8 µg/ml). The optimal conditions should be determined individually. It detects p53 in samples of human, mouse and rat origins. Other applications have not been determined. Blocking peptide (Cat.# PK-AB577-3036P) is available separately.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.