

# p53AIP1 antibody (pAb)

## Rabbit Anti-Human p53AIP1

### Instruction Manual

Catalog Number	PK-AB718-2449
Synonyms	p53AIP1 Antibody; p53AIP1, p53-regulated apoptosis-inducing protein 1, p53AIP
Description	The p53 tumor-suppressor protein can induce apoptosis through transcriptional activation of several genes. One such protein p53AIP was initially identified through direct cloning of p53 binding sequences from human genomic DNA. Its expression is inducible by p53 following p53 phosphorylation on Ser-46, and ectopic expression of p53AIP leads to apoptotic cell death. Both the phosphorylation of p53 and the induction of p53AIP were blocked by inhibiting the expression of p53DINP1 by the introduction of antisense oligonucleotides to p53DINP1, suggesting that the apoptosis associated with p53AIP expression is regulated by p53DINP1. Finally, as adenovirus-mediated introduction of p53AIP has been shown to suppress tumor growth in vivo, it has been suggested that p53AIP gene transfer may become a useful strategy for the treatment of p53-resistant cancers. Three isoforms of p53AIP are known to exist; this antibody will detect all three.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal p53AIP antibody was raised against a 15 amino acid peptide from near the amino terminus of human p53AIP (GenBank accession no. BAB16421).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, ICC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually.  p53AIP1 antibody can be used for detection of p53AIP1 by Western blot at 4 - 8 µg/mL. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Kern SE, Pietenpol JA, Thiagalingam S, Seymour A, et al. Oncogenic forms of p53 inhibit p53-regulated gene expression. Science 1991; 252:1708-11 Oda K, Arakawa H, Tanaka T, et al. p53AIP1, a potential mediator of p53-dependent apoptosis, and its regulation by Ser- 46-phosphorylated p53. Cell 2000; 102:849-52. Okamura S, Arakawa H, Tanaka T, et al. p53DINP1, a p53- inducible gene, regulates p53-dependent apoptosis. Mol. Cell 2001; 8:85-94. 4. Yoshida K, Monden M, Nakamura Y, et al. Adenovirusmediated p53AIP1 gene transfer as a new strategy for treatment of p53-resistant tumors. Cancer Sci. 2004; 95:91-7. (07-01D)
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2449P HL60 Cell Lysate, Cat. No. PK-AB718-1209 p53DINP1 Antibody, Cat. No. PK-AB718-3045
FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	