

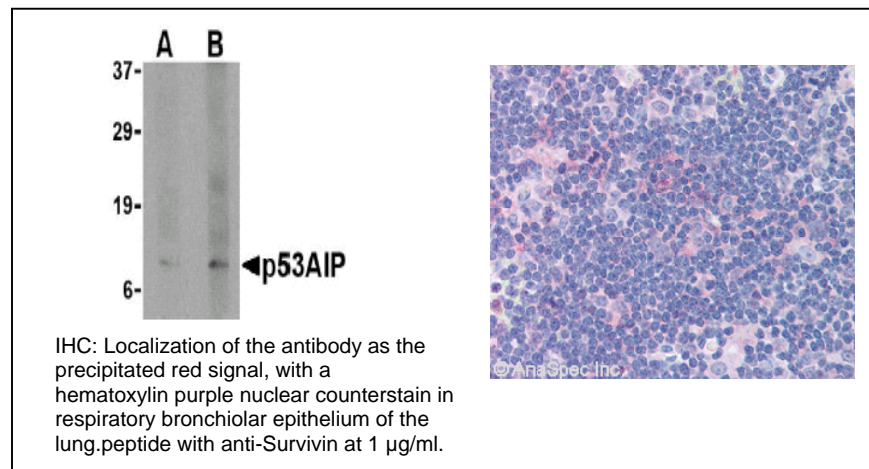


Product Data Sheet

Product Name:	Anti-p53AIP (CT)
Catalog Number:	28013
Lot Number:	See label on vial
Product Description:	This polyclonal antibody is supplied as an epitope affinity purified rabbit IgG, 50 µg in 250 µl (0.2 mg/ml) of 1X PBS (pH 7.4) containing 0.05% sodium azide.
Immunogen:	A synthetic peptide corresponding to C-terminus of human p53AIP1.
Species Reactivity:	This antibody can be used for detection of p53AIP1 by Western blot. An immune reactive band was observed from apoptotic A431 or HEK293 cell lysate. In IHC, it showed moderate to strong staining of a number of cell types in formalin-fixed human tissues including heart, kidney, lung, and pancreas.
Application Notes:	The following concentration ranges are recommended starting points for this product. Optimal working concentrations should be determined by the investigator for specific applications.

Western Blot:	0.5 to 2 µg/ml
Immunohistochemistry :	2.5-20.0 µg/ml (Optimal 20 µg/ml)
IP*:	3.0-5.0 µg/extract from 10 ⁷ cells

(*Recommended but not tested)



Background:

The p53 tumor-suppressor protein can induce apoptosis through transcriptional activation of several genes (1). One such protein p53AIP was initially identified through direct cloning of p53 binding sequences from human genomic DNA (2). 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 (3). 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 (4). Three isoforms of p53AIP are known to exist; this antibody will detect all three.

Storage:

Store at 2-8°C for up to one year. Avoid repeated freezing and thawing.

References:

1. Oda, K. et al. *Cell* **102**, 849 (2000).
2. Okamura, S. et al. *Mol Cell* **8**, 85 (2001).
3. Costanzo, A. et al. *Mol Cell* **9**, 175 (2002).
4. Yamaguchi, T. et al. *Cancer Res* **61**, 8256 (2001).

This product is for *in vitro* research use only.