

p53DINP1 / SIP antibody (pAb)

Rabbit Anti-Human/Mouse/Rat p53DINP1 (Apoptosis Associated p53 Target)

Instruction Manual

Catalog Number	PK-AB718-3045
Synonyms	p53DINP1 Antibody: p53DINP1, SIP
Description	Apoptosis is related to many diseases and development. The p53 tumor-suppressor protein induces apoptosis through transcriptional activation of several genes. A novel p53 inducible gene was identified recently and designated p53DINP1 (for p53-dependent damage-inducible nuclear protein 1) and SIP (for stress induced protein) in human and mouse. A p53DINP1 antisense oligonucleotide inhibits and overexpression of p53DINP1 enhances Ser46 phosphorylation of p53, induction of p53AIP1, and cell death induced by DNA double-strand breaks. p53DINP1 may regulate p53-dependent apoptosis through phosphorylation at Ser46 and induction of p53AIP1. The p53DINP1/SIP gene encodes two proteins of 27 and 18 kDa in human and mouse termed p53DINP1-a and p53DINP1-b or SIP27 and SIP18. p53DINP1/SIP is expressed in many tissues and induced by a variety of stress agents including UV stress, mutagenic stress, heat shock, and oxidative stress.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal p53DINP1 antibody was raised with a synthetic peptide corresponding to 14 amino acids near the amino terminus of human p53DINP1 (GenBank Accession number Q96A56). This sequence is identical between α and β forms of the p53DINP1 proteins, and differs by one amino acid from those of mouse.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	A lower band at 18 kDa was detected in human spleen, and mouse liver and kidney tissue lysates, which may represent the p53DINP1-b form.
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, IHC Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. p53DINP1 antibody can be used for detection of p53DINP1 by Western blot at 0.5 - 1 µg/mL. A band at 27 kDa can be detected. Antibody can also be used for immunohistochemistry starting at 2 µg/mL.
Images	Available upon request.
References	Okamura S, Arakawa H, Tanaka T, Nakanishi H, Ng CC, Taya Y, Monden M, Nakamura Y. p53DINP1, a p53-inducible gene, regulates p53-dependent apoptosis. Mol Cell. 2001;8(1):85-94. Tomasini R, Samir AA, Vaccaro MI, Pebusque MJ, Dagorn JC, Iovanna JL, Dusetti NJ. Molecular and functional characterization of the stress-induced protein (SIP) gene and its two transcripts generated by alternative splicing. SIP induced by stress and promotes cell death. J Biol Chem. 2001;276(47):44185-92.
Images	Available upon request.
Related Products	IN VITRO RES Blocking Peptide, Cat. No. PK-AB718-3045P, Human Tissue Lysate, Cat. No. PK-AB718-1302