

p53R2 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat p53R2 (Novel Ribonucleotide Reductase)

Instruction Manual

| | |
|---------------------|---|
| Catalog Number | PK-AB718-2383 |
| Synonyms | p53R2 Antibody: |
| Description | The p53 tumor-suppresspr gene integrates numerous signals that control cell life and death. Several novel molecules involved in p53 signaling, including p53R2 , Chk2, p53AIP1, Noxa, PIDD, and PID/MTA2, were recently discovered. p53R2 is a p53 inducible gene that contains a p53 binding sequence and encodes a subunit of the enzyme ribonucleotide reductase. p53R2 is induced by the reagents, ultraviolet and γ -irradiation that cause DNA damages. The product of p53R2 gene is directly involved in the p53 checkpoint for repair of damaged DNA. The isoform of the p53 family member p73 also induces p53R2 expression. p53R2 is an important target of p53 for tumour suppression. |
| Quantity | 100 μ g |
| Source / Host | Rabbit |
| Immunogen | Rabbit polyclonal p53R2 antibody was raised against a synthetic peptide (GDPERPEAAGLDQDER) corresponding to amino acids 2 to 17 of human p53R2 . |
| Purification Method | Affinity chromatography purified via peptide column. |
| Clone / IgG Subtype | Polyclonal antibody |
| Species Reactivity | Human, Mouse, Rat |
| 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 μ l 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, 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. p53R2 antibody can be used for detection of p53R2 by Western blot at 0.5 to 1 μ g/mL. A 39 kDa band can be detected. Antibody can also be used for immunohistochemistry starting at 1 μ g/mL. For immunofluorescence start at 20 μ g/mL. |
| Images | Available upon request. |
| References | Matsuoka S, Huang M, Elledge SJ. Linkage of ATM to cell cycle regulation by the Chk2 protein kinase. <i>Science</i> . 1998;282:1893-7. Tanaka H, Arakawa H, Yamaguchi T, Shiraishi K, Fukuda S, Matsui K, Takei Y, Nakamura Y. A ribonucleotide reductase gene involved in a p53-dependent cell-cycle checkpoint for DNA damage. <i>Nature</i> . 2000;404:42-9. Oda E, Ohki R, Murasawa H, Nemoto J, Shibue T, Yamashita T, Tokino T, Taniguchi T, Tanaka N. Noxa, a BH3-only member of the Bcl-2 family and candidate mediator of p53-induced apoptosis. <i>Science</i> . 2000;288(5468):1053-8. Oda K, Arakawa H, Tanaka T, Matsuda K, Tanikawa C, Mori T, Nishimori H, Tamai K, Tokino T, Nakamura Y, Taya Y. p53AIP1, a potential mediator of p53-dependent apoptosis, and its regulation by Ser-46-phosphorylated p53. <i>Cell</i> . 2000 Sep 15;102(6):849-62. |
| Images | Available upon request. |
| Related Products | A431 Lysate, Cat. No. PK-AB718-1202 p53R2 Peptide, Cat. No. PK-AB718-2383P |

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