

MTBP (IN2) antibody (pAb)

Rabbit Anti-Human MTBP (MDM2 Binding Protein)

Instruction Manual

Catalog Number	PK-AB718-2447
Synonyms	MTBP Antibody; MTBP
Description	The p53 tumor-suppressor gene integrates numerous signals that control cell life and death. Several novel molecules involved in p53 network, including Chk2 , p53R2, p53AIP1, Noxa, PIDD, PID/MTA2 and MTBP, were recently discovered. The transcriptional activity of p53 is modulated by posttranslational regulations of the p53 protein including stabilization and acetylation. P53 transcriptionally activates MDM2 gene then the translated MDM2 protein binds to p53 and promotes the degradation of p53 leading to lowering the concentration of p53 protein. MDM2 inhibits both p53 mediated G1 arrest and apoptosis. A recently discovered protein termed MTBP was found to bind to MDM2 and to inhibit the modulation effect of MDM2 on p53. MTBP is expressed in a variety of normal tissues.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal MTBP antibody was raised against a synthetic peptide (GAVECFEEEDSNSRESLS) corresponding to amino acids 122 to 139 of human MTBP, which differs from the mouse sequence by three amino acids .
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. MTBP antibody can be used for detection of MTBP by Western blot at 1 µg/mL. A 104 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/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	K562 Lysate, Cat. No. PK-AB718-1204 MTBP (IN2) Peptide, Cat. No. PK-AB718-2447P

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