

# PID / MTA2 antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat PID (Involved in p53 Deacetylation)

### Instruction Manual

Catalog Number	PK-AB718-2443
Synonyms	PID Antibody: PID, MTA2
Description	The p53 tumor-suppressor gene integrates numerous signals that control cell life and death. Several novel molecules involved in p53 pathway, including Chk2, p53R2, p53AIP1, Noxa, PIDD, and PID/MTA2, were recently discovered. The transcriptional activity of p53 is modulated by protein stability and acetylation. PID/MTA2, also termed MTA1-L1, was found to be a subunit of nucleosome remodeling and deacetylating (NRD/NuRD) complex. PID/MTA2 modulates the enzymatic activity of the histone deacetylase complex and its expression reduces the levels of acetylated p53. Deacetylation of p53 by PID/MTA2 represses p53-dependent transcriptional activation and modulates p53-mediated cell growth arrest and apoptosis. PID/MTA2 is ubiquitously expressed in human tissues.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal PID antibody was raised against a synthetic peptide (PAPSHPASTNEPIVLED) corresponding to amino acids 652 to 668 of human PID/MTA2, which differ from the mouse sequence by one amino acid.
Purification Method	Antibody is purified via DEAE-column chromatography.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	No cross response to MTA1.
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.  PID antibody can be used for detection of PID by Western blot at 1 µg/mL. A 75 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 10 µg/mL. For immunofluorescence start at 10 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	HeLa Lysate, Cat. No. PK-AB718-1201 PID/MTA2 Peptide, Cat. No. PK-AB718-2443P

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