

PERP antibody (pAb)

Rabbit Anti-Human PERP (Mediator of p53 Induced Apoptosis)

Instruction Manual

Catalog Number	PK-AB718-2451
Synonyms	PERP Antibody: PERP
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, MTBP and PERP, were identified and their genes were cloned recently. PERP, also termed PIGPC1 and THW, is a plasma membrane protein. p53 binds to the promoter of PERP and transcriptionally activates PERP gene then the translated PERP protein mediates the p53 induced apoptosis. The expression of PERP causes cell death. PERP is a mediator of p53 induced apoptosis. PERP has sequence similarity to PMP-22/gas3 and is a new member of the PMP-22/gas3 family.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal PERP antibody was raised against a synthetic peptide corresponding to amino acids near the carboxy terminus of human PERP, which differ from the mouse sequence by three amino acids (7-9).
Purification Method	Antibody is purified via DEAE-column chromatography.
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 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. PERP antibody can be used for detection of PERP by Western blot at 1 µg/mL. A 21 kDa band can be detected. Antibody can also be used for immunocytochemistry starting 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	Blocking Peptide, Cat. No. PK-AB718-2451P A431 Cell Lysate, Cat. No. PK-AB718-1202 CHK2 Antibody (NT), Cat. No. PK-AB718-2391 p53R2 Antibody (NT), Cat. No. PK-AB718-2385

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