

# Chk2 (NT) antibody (pAb)

## Rabbit Anti-Human/Mouse/Rat Chk2 (Novel Check Point Kinase)

### Instruction Manual

Catalog Number	PK-AB718-2391
Synonyms	Chk2 Antibody:
Description	The p53 tumor-suppressor gene integrates numerous signals that control cell life and death. Several novel molecules involved in p53 signaling, including Chk2, p53R2, p53AIP1, Noxa, PIDD, and PID/MTA2, were recently discovered. The checkpoint kinase Chk2 is the mammalian homologue of yeast Cds1/Rad53. In response to DNA damage, the checkpoint kinase ATM phosphorylates and activates Chk2, which in turn directly phosphorylates and activates p53. Chk2 serves as ATM downstream effector to mediate activation of p53. Chk2 also phosphorylates and activates BRCA1, the product of a tumor suppressor gene that is mutated in breast and ovarian cancer.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Chk2 antibody was raised against a synthetic peptide corresponding to amino acids near the amino terminus of human Chk2.
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 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. Application have to be determined individually. Chk2 antibody can be used for detection of Chk2 by Western blot at 1 µg/mL. A 60 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 1 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	Chk2 Peptide, Cat. No. PK-AB718-2391P Jurkat Lysate, Cat. No. PK-AB718-1205 Noxa Antibody, Cat. No. PK-AB718-2437

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