

### Instruction Manual

<b>Catalog Number</b>	PK-AB718-2437
<b>Synonyms</b>	Noxa Antibody: Noxa, PMA-induced protein 1, PMAIP1, APR
<b>Description</b>	Apoptosis is related to many diseases and development. The p53 tumor-suppressor protein induces apoptosis through transcriptional activation of several genes including p53R2, p53AIP1, and PUMA. A new p53 target gene, Noxa, was recently identified, which encodes a protein belonging to the subfamily of BH3-only proapoptotic proteins. Noxa and PUMA are both transcriptional targets of p53 and BH3-only proteins. X-ray irradiation increased p53-dependent Noxa mRNA and protein levels. Noxa, when ectopically expressed, interacted with anti-apoptotic Bcl-2 family members, resulting in the activation of caspase-9. Noxa, like PUMA, localized to mitochondria and induces apoptosis in response to p53. Noxa and PUMA may represent direct mediators of p53-induced apoptosis. Increased levels of p53 and its target gene Noxa was found in the impaired tumor development.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	Rabbit Noxa polyclonal antibody was raised against a synthetic peptide corresponding to 17 amino acids at the amino terminus of mouse Noxa (GenBank accession no. NP_067426).
<b>Purification Method</b>	Immunoaffinity chromatography purified IgG.
<b>Clone / IgG Subtype</b>	Polyclonal antibody
<b>Species Reactivity</b>	Human, Mouse, Rat
<b>Specificity</b>	
<b>Formulation</b>	Antibody is supplied in PBS containing 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	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.
<b>Applications</b>	E, WB, IHC, IF <span style="float: right;">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.</span>  Noxa antibody can be used for detection of Noxa by Western blot at at 0.5 - 2 µg/mL. Antibody can also be used for immunohistochemistry starting at 1 µg/mL. For immunofluorescence start at 10 µg/mL.
<b>Images</b>	Available upon request.
<b>References</b>	Oda E, Ohki R, Murasawa H, et al. Noxa, a BH3-only member of the Bcl-2 family and candidate mediator of p53-induced apoptosis. <i>Science</i> 2000; 288:1053-8. Nakano K and Vousden KH. PUMA, a novel proapoptotic gene, is induced by p53. <i>Mol. Cell</i> 2001; 7:683-94. Yu J, Zhang L, Hwang PM, et al. PUMA induces the rapid apoptosis of colorectal cancer cells. <i>Mol. Cell</i> 2001; 7:673-82. Eferl R, Ricci R, Kenner L, et al. Liver tumor development. c-Jun antagonizes the proapoptotic activity of p53. <i>Cell</i> 2003; 112:181-92.
<b>Images</b>	Available upon request.
<b>Related Products</b>	Stomach Lysate, Catalog No. PK-AB718-1344-N; Noxa Peptide, Cat. No. PK-AB718-2437P Bmf Antibody, Cat. No. PK-AB718-3031; Bim Antibody, Cat. No. PK-AB718-2065 p53DINP1 Antibody, Cat. No. PK-AB718-3045; Chk2 Antibody, Cat. No. PK-AB718-2391; p53R2 Antibody, Cat. No. PK-AB718-2383; PUMA Antibody, Cat. No. PK-AB718-3041
FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.	