

# NIK (CT) antibody (pAb)

## Rabbit Anti-Human NF-κB Inducing Kinase

### Instruction Manual

|                     |   |
|---------------------|---|
| Catalog Number      | PK-AB718-1129   |
| Synonyms            | NIK Antibody:   |
| Description         | Nuclear factor κ B (NF-κB) is a ubiquitous transcription factor and an essential mediator of gene expression during activation of immune and inflammatory responses. NF-κB mediates the expression of a great variety of genes in response to extracellular stimuli including IL-1, TNFα, LPS and mitogens. A serine/threonine protein kinase which mediates NF-κB activation by IL-1, TNFα and CD95 was identified recently and designated NIK (for NF-κB inducing kinase) . NIK is an activator of IκB kinase α and β (IKKα and IKKβ). Therefore, NIK is a key molecule in the NF-κB signaling pathway leading to the induction of a variety of gene expression in response to proinflammatory cytokines and bacteria products. |
| Quantity            | 100 µg  |
| Source / Host       | Rabbit  |
| Immunogen           | Rabbit polyclonal NIK antibody was raised against a peptide corresponding to amino acids 931 to 947 of human NIK .  |
| 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<br>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.   |
|                     | NIK antibody can be used for detection of NIK by Western blot.  |
| Images              | NA  |
| References          | Malinin NL, Boldin MP, Kovalenko AV, Wallach . MAP3K-related kinase involved in NF-κB induction by TNF, CD95 and IL-1. Nature 1997;385:540-544<br>Regnier CH, Song HY, Gao X, Goeddel DV, Cao Z, Rothe M. Identification and characterization of an IκB kinase. Cell 1997;90:373-383<br>Woronicz JD, Gao X, Cao Z, Rothe M, Goeddel DV. IκB kinase-β: NF-κB activation and complex formation with IκB kinase-α and NIK. Science 1997;278:866-869<br>Ling L, Cao Z, Goeddel D. NF-κB-inducing kinase activates IKK-α by phosphorylation of Ser-176. Proc Natl Acad Sci USA 1998;95:3792-3797   |
| Images              | NA  |
| Related Products    | 293 Lysate, Cat. No. PK-AB718-1209<br>NIK Peptide, Cat. No. PK-AB718-1129P  |

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