

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-2025P
<b>Quantity</b>	50 µg
<b>Source</b>	19 amino acids near the carboxy terminus of human IKK alpha
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 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>	Store IKK alpha peptide at -20°C, stable for one year.
<b>Application</b>	IKK alpha peptide is used for blocking the activity of IKK alpha antibody.
<b>References</b>	<p>DiDonato JA, Hayakawa M, Rothwarf DM, Zandi E, Karin M. A cytokine-responsive IκappaB kinase that activates the transcription factor NF-kappaB. <i>Nature</i> 1997;388:548-54.</p> <p>Liu X, Li P, Widlak P, Zou H, Luo X, Garrard WT, Wang X The 40-kDa subunit of DNA fragmentation factor induces DNA fragmentation and chromatin condensation during apoptosis. <i>Proc Natl Acad Sci USA</i> 1998;95:8461-6</p> <p>Halenbeck R, MacDonald H, Roulston A, Chen TT, Conroy L, Williams LT. CPAN, a human nuclease regulated by the caspase-sensitive inhibitor DFF45. <i>Curr Biol</i> 1998;8:537-40</p>

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