

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-2329P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the carboxy terminus of human IKK epsilon
<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 epsilon peptide at -20°C, stable for one year.
<b>Application</b>	IKK epsilon peptide is used for blocking the activity of IKK epsilon antibody.
<b>References</b>	<p>Peters RT, Liao SM, Maniatis T. IKKepsilon is part of a novel PMA-inducible IκappaB kinase complex. <i>Mol Cell</i> 2000;5(3):513-22.</p> <p>Sheridan JP, Marsters SA, Pitti RM, et al. A. Control of TRAIL-induced apoptosis by a family of signaling and decoy receptors. <i>Science</i> 1997;277:818-21</p> <p>Degli-Esposti MA, Smolak PJ, Walczak H, et al, Smith CA. Cloning and characterization of TRAIL-R3, a novel member of the emerging TRAIL receptor family. <i>J Exp Med</i> 1997;186(7):1165-70</p>

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