

Instruction Manual

Catalog Number	PK-AB718-2121P
Quantity	50 µg
Source	19 amino acids near the carboxy terminus of human IKK beta (C3)
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 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	Store IKK beta (C3) peptide at -20°C, stable for one year.
Application	IKK beta peptide is used for blocking the activity of IKK beta antibody.
References	<p>Zandi E, Rothwarf DM, Delhase M, Hayakawa M, Karin M. The IκappaB kinase complex (IKK) contains two kinase subunits, IKKα and IKKβ, necessary for IκappaB phosphorylation and NF-κappaB activation. <i>Cell</i> 1997;91:243-52.</p> <p>Samson M, Stordeur P, Labbe O, Soularue P, Vassart G, Parmentier M. Molecular cloning and chromosomal mapping of a novel human gene, ChemR1, expressed in T lymphocytes and polymorphonuclear cells and encoding a putative chemokine receptor. <i>Eur J Immunol</i> 1996;26:3021-8</p> <p>Goya I, Gutierrez J, Varona R, Kremer L, Zaballos A, Marquez G. Identification of CCR8 as the specific receptor for the human beta-chemokine I-309: cloning and molecular characterization of murine CCR8 as the receptor for TCA-3. <i>J Immunol</i> 1998;160:1975-81</p>

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