

Instruction Manual

Catalog Number	PK-AB718-2115P
Quantity	50 µg
Source	17 amino acids near the carboxy terminus of human IKK alpha (C2)
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 alpha (C2) peptide at -20°C, stable for one year.
Application	IKK alpha peptide is used for blocking the activity of IKK alpha antibody.
References	DiDonato JA, Hayakawa M, Rothwarf DM, Zandi E, Karin M. A cytokine-responsive IκappaB kinase that activates the transcription factor NF-kappaB. Nature 1997;388:548-54.

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