

Instruction Manual

| | |
|--------------------------------|---|
| Catalog Number | PK-AB718-2335P |
| Quantity | 50 µg |
| Source | 17 amino acids near the carboxy terminus of human IKK gamma |
| 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 gamma peptide at -20°C, stable for one year. |
| Application | IKK gamma peptide is used for blocking the activity of IKK gamma antibody. |
| References | Rothwarf DM, Zandi E, Natoli G, Karin M. IKK-gamma is an essential regulatory subunit of the IκappaB kinase complex. Nature 1998;395(6699):297-300. |

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