

Instruction Manual

Catalog Number	PK-AB718-3395P
Quantity	50 µg
Source	16 amino acids near the amino terminus of human BTK (NT)
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 BTK peptide at -20°C, stable for one year.
Application	Casp-6 peptide is used for blocking the activity of BTK antibody.
References	<p>Vetrie D, Vorechovsky I, Sideras P, et al. The gene involved in X-linked agammaglobulinaemia is a member of the src family of tyrosine kinases. <i>Nature</i> 1993; 361:226-33.</p> <p>Deveraux QL, Leo E, Stennicke HR, et al. Cleavage of human inhibitor of apoptosis protein XIAP results in fragments with distinct specificities for caspases. <i>EMBO J.</i> 1999; 18:5242-51.</p> <p>Holcik M, Yeh C, Korneluk RG, et al. Translational upregulation of X-linked inhibitor of apoptosis (XIAP) increases resistance to radiation induced cell death. <i>Oncogene</i> 2000; 19:4174-7.</p> <p>Hockenbery D, Nunez G, Milliman C, et al. Bcl-2 is an inner mitochondrial membrane protein that blocks programmed cell death. <i>Nature</i> 1990; 348:334-6.</p>

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