

Instruction Manual

Catalog Number	PK-AB718-3355P
Quantity	50 µg
Source	14 amino acids near the center of human Bid
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 Bid peptide at -20°C, stable for one year.
Application	Bid peptide is used for blocking the activity of Bid antibody.
References	<p>Wang K, Yin XM, Chao DT, et al. BID: a novel BH3 domain-only death agonist. <i>Genes Dev.</i> 1996; 10:2859-69.</p> <p>Vilas GL, Corvi MM, Plummer GJ, et al. Posttranslational myristoylation of caspase-activated p21-activated protein kinase 2 (PAK2) potentiates late apoptotic events. <i>Proc. Natl. Acad. Sci. USA</i> 2006; 103:6542-7.</p> <p>Janeway CA Jr and Medzhitov R. Innate immune recognition. <i>Annu. Rev. Immunol.</i> 2002; 20:197-216.</p>

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