

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3819P
<b>Quantity</b>	50 µg
<b>Source</b>	15 amino acids near the amino terminus of human Bik
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	Store Bik peptide at -20°C, stable for one year.
<b>Application</b>	Bik peptide is used for blocking the activity of Bik antibody.
<b>References</b>	<p>Germain M, Mathai JP, and Shore GC. BH-3-only BIK functions at the endoplasmic reticulum to stimulate cytochrome c release from mitochondria. <i>J. Biol. Chem.</i> 277:18053-60.</p> <p>Haruta I, Kato Y, Hashimoto E, et al. Association of AIM, a novel apoptosis inhibitory factor, with hepatitis via supporting macrophage survival and enhancing phagocytic function of macrophages. <i>J. Biol. Chem.</i> 2001; 276:22910-4</p> <p>Kuwata K, Watanabe H, Jiang S-Y, et al. AIM inhibits apoptosis of T cells and NKT cells in Corynebacterium-induced granuloma formation in mice. <i>Am. J. Path.</i> 2003; 162:837-47.</p> <p>Meyers JH, Chakravarti S, Schlesinger D, et al. TIM-4 is the ligand for TIM-1, and the TIM-1-TIM4 interaction regulates T cell proliferation. <i>Nat. Immunol.</i> 2005; 6:455-64.</p>

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