

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3351P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the amino terminus of human Bax
<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 Bax peptide at -20°C, stable for one year.
<b>Application</b>	Bax peptide is used for blocking the activity of Bax antibody.
<b>References</b>	Oltvai ZN, Milliman CL, and Korsmeyer SJ. Bcl-2 heterodimerizes in vivo with a conserved homolog, Bax, that accelerates programmed cell death. <i>Cell</i> 1993; 74:609-19. Janeway CA Jr. and Medzhitov R. Innate immune recognition. <i>Annu. Rev. Immunol.</i> 2002; 20:197-216. O'Neill LAJ, Fitzgerald FA, and Bowie AG. The Toll-IL-1 receptor adaptor family grows to five members. <i>Trends in Imm.</i> 2003; 24:286-9.

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