

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-2065P
<b>Quantity</b>	50 µg
<b>Source</b>	19 amino acids near the center of human Bim
<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 Bim peptide at -20°C, stable for one year.
<b>Application</b>	Bim peptide is used for blocking the activity of Bim antibody.
<b>References</b>	<p>O'Connor,L., Strasser,A., O'Reilly,L.A., Hausmann,G., Adams,J.M., Cory,S. and Huang,D.C. Bim: a novel member of the Bcl-2 family that promotes apoptosis. <i>EMBO J.</i> 1998;17:384-395.</p> <p>Sheridan JP, Marsters SA, Pitti RM, Gurney A, Skubatch M, Baldwin D, Ramakrishnan L, Gray CL, Baker K, Wood WI, Goddard AD, Godowski P, Ashkenazi A. Control of TRAIL-induced apoptosis by a family of signaling and decoy receptors. <i>Science</i> 1997;277:818-21</p> <p>Marsters SA, Sheridan JP, Pitti RM, Huang A, Skubatch M, Baldwin D, Yuan J, Gurney A, Goddard AD, Godowski P, Ashkenazi A. A novel receptor for Apo2L/TRAIL contains a truncated death domain. <i>Curr Biol</i> 1997;7:1003-6</p>

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