

Instruction Manual

Catalog Number	PK-AB718-3405P
Quantity	50 µg
Source	16 amino acids near the middle of human Bim
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 Bim peptide at -20°C, stable for one year.
Application	Bim peptide is used for blocking the activity of Bim antibody.
References	<p>O'Connor L, Strasser A, O'Reilly LA, et al. Bim: a novel member of the Bcl-2 family that promotes apoptosis. <i>EMBO J.</i> 1998; 17:384-395.</p> <p>Ottile S, Diaz JL, Horne W, et al. Dimerization properties of human BAD. Identification of a BH-3 domain and analysis of its binding to mutant BCL-2 and BCL-XL proteins. <i>J. Biol. Chem.</i> 1997; 272:30866-72.</p> <p>Zhou XM, Liu Y, Payne G, et al. Growth factors inactivate the cell death promoter BAD by phosphorylation of its BH3 domain on Ser155. <i>J. Biol. Chem.</i> 2000; 275:25046-51.</p> <p>Kiefer M, Brauer MJ, Powers VC, et al. Modulation of apoptosis by the widely distributed Bcl-2 homologue Bak. <i>Nature</i> 1995; 374:736-9.</p>

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