

Instruction Manual

Catalog Number	PK-AB718-3165P
Quantity	50 µg
Source	15 amino acids near the amino terminus of human Bcl-G
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 Bcl-G peptide at -20°C, stable for one year.
Application	Bcl-G peptide is used for blocking the activity of Anit-Bcl-G antibody.
References	<p>Guo B, Godzik A, and Reed JC. Bcl-G, a novel pro-apoptotic member of the Bcl-2 family. <i>J. Biol. Chem.</i> 2000; 276:2780-5.</p> <p>Shu HB, Hu WH, Johnson H. TALL-1 is a novel member of the TNF family that is down-regulated by mitogens. <i>J Leukoc Biol</i> 1999;65:680-3</p> <p>Mukhopadhyay A, Ni J, Zhai Y, Yu GL, Aggarwal BB. Identification and characterization of a novel cytokine, THANK, a TNF homologue that activates apoptosis, nuclear factor-kappaB, and c-Jun NH2-terminal kinase. <i>J Biol Chem</i> 1999 ;274:15978-81</p>

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