

Instruction Manual

Catalog Number	PK-AB718-3337P
Quantity	50 µg
Source	14 amino acids near the center of human Bcl-2
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-2 peptide at -20°C, stable for one year.
Application	Bcl-2 peptide is used for blocking the activity of Bcl-2 antibody.
References	<p>Hockenbery D, Nunez G, Milliman C, et al. Bcl-2 is an inner mitochondrial membrane protein that blocks programmed cell death. <i>Nature</i> 1990; 348:334-6.</p> <p>Bieniasz PD. Intrinsic immunity: a front-line defense against viral attack. <i>Nat Immunol.</i> 2004; 5:1109-15.</p> <p>Sheehy AM, Gaddis NC, Choi JD, et al. Isolation of a human gene that inhibits HIV-1 infection and is suppressed by the viral Vif protein. <i>Nature</i> 2002; 418:646-50.</p> <p>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.</p>

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