

Instruction Manual

Catalog Number	PK-AB718-3023P
Quantity	50 µg
Source	15 amino acids near the center of human Bcl-rambo
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-rambo peptide at -20°C, stable for one year.
Application	Bcl-rambo peptide is used for blocking the activity of the Bcl-rambo antibody.
References	<p>Kataoka T, Holler N, Michau O, et al. Bcl-rambo, a novel Bcl-2 homologue that induces apoptosis via its unique C-terminal extension. <i>J. Biol. Chem.</i> 2001; 276:19548-54.</p> <p>Hugot JP, Chamaillard M, et al. Association of NOD2 leucine-rich repeat variants with susceptibility to Crohn's disease. <i>Nature</i> 2001; 411:599-603.</p> <p>Ogura Y, Bonen DK, Inohara N, et al. A frameshift mutation in NOD2 associated with susceptibility to Crohn's disease. <i>Nature</i> 2001; 411:603-6.</p>

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