

Instruction Manual

Catalog Number	PK-AB718-3873P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human Bfl-1
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 Bfl-1 peptide at -20°C, stable for one year.
Application	Bfl-1 peptide is used for blocking the activity of Bfl-1 antibody.
References	<p>Choi SS, Park IC, Yun JW, et al. A novel Bcl-2 related gene, Bfl-1, is overexpressed in stomach cancer and preferentially expressed in bone marrow. <i>Oncogene</i> 1995; 11:1693-8.</p> <p>Sinha SK and Chaudhary PM. Induction of apoptosis by X-linked ectodermal dysplasia receptor via a caspase 8-dependent mechanism. <i>J. Biol. Chem.</i> 2004; 41873-81.</p> <p>Martinon F and Tschopp J. NLRs join TLRs as innate sensors of pathogens. <i>TRENDS Imm.</i> 2005; 26:447-54.</p>

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