

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3991P
<b>Quantity</b>	50 µg
<b>Source</b>	17 amino acids near the amino terminus of human Blimp-1
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	Store Blimp-1 peptide at -20°C, stable for one year.
<b>Application</b>	Blimp-1 peptide is used for blocking the activity of Blimp-1 antibody.
<b>References</b>	Turner CAJ, Mack DH, and Davis MM. Blimp-1, a novel zinc finger-containing protein that can drive the maturation of B-lymphocytes into immunoglobulin-secreting cells. Cell 1994; 77:297-306.

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