

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3097P
<b>Quantity</b>	50 µg
<b>Source</b>	15 amino acids near the carboxy terminus of human BAFF Receptor
<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 BAFF Receptor peptide at -20°C, stable for one year.
<b>Application</b>	BAFF receptor peptide is used for blocking the activity of BAFF-R antibody.
<b>References</b>	Thompson JS, Bixler SA, Qian F, et al. BAFF-R, a newly identified TNF receptor that specifically interacts with BAFF. <i>Science</i> . 2001;293:2108-11. Agostini L, Martinon F, Burns K, et al. NALP3 forms an IL-1beta-processing inflammasome with increased activity in Muckle-Wells autoinflammatory disorder. <i>Immunity</i> 2004; 20:319-25. Martinon F and Tschopp J. NLRs join TLRs as innate sensors of pathogens. <i>TRENDS Imm.</i> 2005; 26:447-54.

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