

Instruction Manual

Catalog Number	PK-AB718-2221P
Quantity	50 µg
Source	16 amino acids near the carboxy terminus of human BAFF
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 BAFF peptide at -20°C, stable for one year.
Application	BAFF peptide is used for blocking the activity of BAFF antibody.
References	<p>Moore PA, Belvedere O, Orr A, et al. BLYS: member of the tumor necrosis factor family and B lymphocyte stimulator. <i>Science</i> 1999;285:260-3.</p> <p>Alkhatib G, Liao F, Berger EA, Farber JM, Peden KW. A new SIV co-receptor, STRL33. <i>Nature</i> 1997;388:238</p> <p>Loetscher M, Amara A, Oberlin E, Brass N, Legler D, Loetscher P, D'Apuzzo M, Meese E, Rousset D, Virelizier JL, Baggiolini M, Arenzana-Seisdedos F, Moser B. TYMSTR, a putative chemokine receptor selectively expressed in activated T cells, exhibits HIV-1 coreceptor function. <i>Curr Biol</i> 1997;7:652-60 (RD0400)</p>

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