

Instruction Manual

Catalog Number	PK-AB718-3817P
Quantity	50 µg
Source	15 amino acids near the carboxy terminus of human Bif
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 Bif peptide at -20°C, stable for one year.
Application	Bif peptide is used for blocking the activity of Bif antibody.
References	<p>Cuddeback SM, Yamaguchi H, Komatsu K, et al. Molecular cloning and characterization of bif-1. <i>J. Biol. Chem.</i> 2001; 276:20559-65.</p> <p>Haruta I, Kato Y, Hashimoto E, et al. Association of AIM, a novel apoptosis inhibitory factor, with hepatitis via supporting macrophage survival and enhancing phagocytic function of macrophages. <i>J. Biol. Chem.</i> 2001; 276:22910-</p> <p>Kuwata K, Watanabe H, Jiang S-Y, et al. AIM inhibits apoptosis of T cells and NKT cells in Corynebacterium-induced granuloma formation in mice. <i>Am. J. Path.</i> 2003; 162:837-47.</p>

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