

Instruction Manual

Catalog Number	PK-AB718-3807P
Quantity	50 µg
Source	16 amino acids near the amino terminus of human AIM
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 AIM peptide at -20°C, stable for one year.
Application	AIM peptide is used for blocking activity of AIM antibody.
References	Biancone L, Bowen MA, Lim A, et al. Identification of a novel inducible cell-surface ligand of CD5 on activated lymphocytes. <i>J. Exp. Med.</i> 1996; 184:811-9.

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