

Instruction Manual

Catalog Number	PK-AB718-4539P
Quantity	50 µg
Source	16 amino acids near the amino terminus of human AIFM3
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 AIFM3 peptide at -20°C, stable for one year.
Application	AIFM3 peptide is used for blocking the activity of AIFM3 antibody.
References	Xie Q, Lin T, Zhang Y, et al. Molecular cloning and characterization of a human AIF-like gene with the ability to induce apoptosis. J. Biol. Chem. 2005; 280:19673-81.

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