

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4541P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the center of human AIFM3
<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 AIFM3 peptide at -20°C, stable for one year.
<b>Application</b>	AIFM3 peptide is used for blocking the activity of AIFM3 antibody.
<b>References</b>	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.