

Instruction Manual

Catalog Number	PK-AB718-2301P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human AIF
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 AIF peptide at -20°C, stable for one year.
Application	AIF peptide is used for blocking the activity of AIF antibody.
References	<p>Susin SA, Lorenzo HK, Zamzami N, et al. Molecular characterization of mitochondrial apoptosis-inducing factor. <i>Nature</i> 1999;397:441-6.</p> <p>Chan SL, Yee KS, Tan KM, Yu VC. The Caenorhabditis elegans sex determination protein FEM-1 is a CED-3 substrate that associates with CED-4 and mediates apoptosis in mammalian cells. <i>J Biol Chem.</i> 2000;275(24):17925-8. (WD0102)</p>

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