

Instruction Manual

Catalog Number	PK-AB718-3039P
Quantity	50 µg
Source	12 amino acids near the carboxy terminus of human AATF
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 AATF peptide at -20°C, stable for one year.
Application	AATF peptide is used for blocking the activity of AATF antibody.
References	<p>Page G, Lodige I, Kogel D et al. AATF, a novel transcription factor that interacts with Dlk/ZIP kinase and interferes with apoptosis. FEBS Lett. 1999; 462:187-191.</p> <p>Kamada Y, Funakoshi T, Shintani T, et al. Tor-mediated induction of autophagy via Apg1 protein kinase complex. J. Cell. Biol. 2000; 150:1507-13.</p> <p>Mizushima N, Noda T, Yoshimori T, et al. A protein conjugation system essential for autophagy. Nature 1998; 395:395-8.</p>

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