

Instruction Manual

Catalog Number	PK-AB718-3607P
Quantity	50 µg
Source	16 amino acids near the carboxy terminus of human AES
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 AES peptide at -20°C, stable for one year.
Application	AES peptide is used for blocking the activity of AES antibody.
References	<p>Miyasaka H, Choudhury BK, Hou WE, et al. Molecular cloning and expression of mouse and human cDNA encoding AES and ESG proteins with strong similarity to Drosophila enhancer of split groucho protein. <i>Eur. J. Biochem.</i> 1993; 216:343-52.</p> <p>Wong GW, Wang J, Hug C, et al. A family of Acrp30/ adiponectin structural and functional paralogs. <i>Proc. Natl. Acad. Sci. USA</i> 2004; 101:10302-7.</p> <p>Shapiro L and Scherer PE. The crystal structure of a complement-1q family protein suggests an evolutionary link to tumor necrosis factor. <i>Curr. Biol.</i> 1998; 8:335-8.</p>

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