

Instruction Manual

Catalog Number	PK-AB718-3617P
Quantity	50 µg
Source	17 amino acids near the amino terminus of human APG7
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 APG7 peptide at -20°C, stable for one year.
Application	APG7 peptide is used for blocking the activity of APG7 antibody.
References	<p>Mizushima N, Noda T, Yoshimori T, et al. A protein conjugation system essential for autophagy. <i>Nature</i> 1998; 395:395-8.</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.