

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3611P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the carboxy terminus of human Beclin-1
<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 Beclin-1 peptide at -20°C, stable for one year.
<b>Application</b>	Beclin-1 peptide is used for blocking the activity of Beclin-1 antibody.
<b>References</b>	<p>Liang XH, Kleeman LK, Jiang HH, et al. Protection against fatal Sindbis virus encephalitis by Beclin, a novel Bcl-2-interacting protein. <i>J. Virol.</i> 1998; 72:8586-96.</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.