

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-4427P
<b>Quantity</b>	50 µg
<b>Source</b>	18 amino acids near the center of human ATG16
<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 ATG16 peptide at -20°C, stable for one year.
<b>Application</b>	ATG16 peptide is used for blocking the activity of ATG16 antibody.
<b>References</b>	Mizushima N, Noda T, and Ohsumi Y. Apg16p is required for the function of the apg12p-apg5p conjugate in the yeast autophagy pathway. EMBO J. 1999; 18:3888-96.

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