

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3495P
<b>Quantity</b>	50 µg
<b>Source</b>	14 amino acids near the carboxy terminus of human GβL
<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 GβL peptide at -20°C, stable for one year.
<b>Application</b>	GβL peptide is used for blocking the activity of GβL antibody.
<b>References</b>	<p>Kim D-H, Sarbassov DD, Ali SM, et al. GβL, a positive regulator of the Rapamycin-sensitive pathway required for the nutrient-sensitive interaction between Raptor and mTOR. <i>Mol. Cell</i> 2003; 11:895-904.</p> <p>Gracie JA, Robertson SE, and McInnes IB. Interleukin-18. <i>J. Leukoc. Biol.</i> 2003; 73:213-224.</p> <p>Kamens J, Paskind M, Hugunin M, et al. Identification and characterization of ICH-2, a novel member of the interleukin-1 beta-converting enzyme family of cysteine proteases. <i>J. Biol. Chem.</i> 1995; 270:15250-6.</p> <p>Martinon F, Burns K, and Tschopp J. The inflammasome: a molecular platform triggering activation of the inflammatory caspases and processing of proIL-1beta. <i>Mol. Cell.</i> 2002; 10:417-26.</p>

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