

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3489P
<b>Quantity</b>	50 µg
<b>Source</b>	15 amino acids near the carboxy terminus of mouse Raptor
<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 Raptor peptide at -20°C, stable for one year.
<b>Application</b>	Raptor peptide is used for blocking the activity of the Raptor antibody.
<b>References</b>	<p>Yonezawa K, Tokunaga C, Oshiro N, et al. Raptor, a binding partner of target of rapamycin. <i>Biochem. Biophys. Res. Commun.</i> 2004; 313:437-441.</p> <p>Chu JJ, Rajamanonmani R, Li J, et al. Inhibition of West Nile virus entry by using a recombinant domain III from the envelope glycoprotein. <i>J. Gen. Virol.</i> 2005; 86:405-12.</p> <p>Chu JJ and Ng ML. Interaction of West Nile virus with alpha v beta 3 integrin mediates virus entry into cells. <i>J. Biol. Chem.</i> 2004; 279:54533-41.</p> <p>Gracie JA, Robertson SE, and McInnes IB. Interleukin-18. <i>J. Leukoc. Biol.</i> 2003; 73:213-224.</p>

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