

Instruction Manual

Catalog Number	PK-AB718-3491P
Quantity	50 µg
Source	13 amino acids near the amino terminus of human Raptor
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 Raptor peptide at -20°C, stable for one year.
Application	Raptor peptide is used for blocking the activity of the Raptor antibody.
References	<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>Kuida K, Lippke JA, Ku G, et al. Altered cytokine export and apoptosis in mice deficient in interleukin-1 beta converting enzyme. <i>Science</i> 1995; 267:2000-3.</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>

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