

Instruction Manual

Catalog Number	PK-AB718-3501P
Quantity	50 µg
Source	15 amino acids near the center of human Rheb
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 Rheb peptide at -20°C, stable for one year.
Application	Rheb peptide is used for blocking the activity of the Rheb antibody.
References	Yamagata K, Sanders LK, Kaufman WE, et al. rheb, a growth factor- and synaptic activity-regulated gene, encodes a novel Ras-related protein. <i>J. Biol. Chem.</i> 1994; 269:16333-9. Gracie JA, Robertson SE, and McInnes IB. Interleukin-18. <i>J. Leukoc. Biol.</i> 2003; 73:213-224. Martinon F, Burns K, and Tschopp J. The inflammasome: a molecular platform triggering activation of the inflammatory caspases and processing of proIL-1β. <i>Mol. Cell.</i> 2002; 10:417-26. Gracie JA, Robertson SE, and McInnes IB. Interleukin-18. <i>J. Leukoc. Biol.</i> 2003; 73:213-224.

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