

Instruction Manual

Catalog Number	PK-AB718-4569P
Quantity	50 µg
Source	17 amino acids near the amino terminus of human RGP1
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 RGP1 peptide at -20°C, stable for one year.
Application	RGP1 peptide is used for blocking the activity of RGP1 antibody.
References	Siniossoglou S, Peak-Chew SY, and Pelham HRB. Ric1p and Rgp1 form a complex that catalyzes nucleotide exchange on Ypt6p. EMBO J. 2000; 19:4885-94.

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