

RGP1 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat RGP1 (Retrograde golgi transport homolog 1)

Instruction Manual

Catalog Number	PK-AB718-4569
Synonyms	RGP1 Antibody: Retrograde golgi transport homolog1
Description	Retrograde golgi transport homolog 1 (RGP1) is the mammalian homolog to the yeast RGP1, a protein that forms a tight complex with RIC1. This complex binds Ypt6p and stimulates guanine nucleotide exchange. RGP1 is localized to the Golgi and is thought to be a potential Golgi recycling factor. Rgp1 yeast mutants exhibit defects in retrograde trafficking similar to those seen in yeast with mutations in other retrograde Golgi transport proteins. It is expected that RGP1 plays a similar role in mammalian cells to that seen in yeast.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	RGP1 antibody was raised in rabbits against a 17 amino acid peptide near the amino terminus of the human RGP1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 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	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. RGP1 antibody can be used for detection of RGP1 by Western blot at 1 - 2 µg/mL.
Images	Available upon request.
References	Siniosoglou 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. Panek HR, Conibear E, Bryan JD, et al. Identification of Rgp1, a novel Golgi recycling factor, as a protein required for efficient localization of yeast casein kinase 1 to the plasma membrane. J. Cell Sci. 113:4545-55.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4569P; RGP1 Peptide Cat. No. PK-AB718-1301; Human Heart Tissue Lysate

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