

RGPD5 (NT) antibody (pAb)

Rabbit Anti-Human RGPD5 (RANBP2-like & GRIP domain-containing protein 5, RGP5)

Instruction Manual

| | |
|---------------------|--|
| Catalog Number | PK-AB718-4585 |
| Synonyms | RGPD5 Antibody; RANBP2-like and GRIP domain-containing protein 5, RGP5, RANBP2L1 |
| Description | The RANBP2-like and GRIP domain containing 5 protein (RGPD5) has high similarity to RANBP2, a large RAN-binding protein localized at the cytoplasmic side of the nuclear pore complex. The gene coding for RGPD5 is thought to have arisen from a gene duplication event of RANBP2 as these highly homologous genes are located close to each other at chromosome 2q11-q12. RGPD5 was identified as an HIV dependency factor (HDF), suggesting that RGPD5 may be an important drug target in HIV treatment. At least two isoforms of RGPD5 are known to exist, of which the shorter isoform is expressed primarily in testis, while the longer of the two is expressed at low levels in a number of somatic tissues. |
| Quantity | 100 µg |
| Source / Host | Rabbit |
| Immunogen | RGPD5 antibody was raised against a 13 amino acid peptide from near the amino terminus of human RGPD5. |
| Purification Method | Affinity chromatography purified via peptide column. |
| Clone / IgG Subtype | Polyclonal antibody |
| Species Reactivity | Human |
| 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, IHC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. RGPD5 antibody can be used for detection of RGPD5 by Western blot at 1 µg/mL. Antibody can also be used for immunohistochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL. |
| Images | Available upon request. |
| References | Nothwang HG, Rensing C, Kubler M, et al. Identification of a novel Ran binding protein 2 related gene (RANBP2L1) and detection of a gene cluster on chromosome 2q11-q12. Genomics 1998; 47:383-92. Brass AL, Dykxhoorn DM, Benita Y, et al. Identification of host proteins required for HIV infection through a functional genomic screen. Science 2008; 319:921-6. Wang LF, Zhu HD, Miao SY, et al. Molecular cloning and characterization of a novel testis-specific nucleoporin-related gene. Arch. Androl. 1999; 42:71-84. |
| Images | Available upon request. |
| Related Products | Cat.No. PK-AB718-4585P; RGPD5 Peptide Cat. No. PK-AB718-1287; EL4 Cell Lysate |

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