

GRTP1 (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat GRTP1 (Growth hormone regulated TBC protein 1)

Instruction Manual

Catalog Number	PK-AB718-4589
Synonyms	GRTP1 Antibody: Growth hormone regulated TBC protein 1, TBC1 domain family member 6
Description	Growth hormone regulated TBC protein 1 (GRTP1) was identified through microarray analysis as a gene whose expression is regulated by growth hormone (GH). It contains the TBC signature motif of GTPase activator proteins of Rab-like small GTPases. GRTP1 mRNA is expressed at highest levels in testes, and is also present in kidney, lung, liver, and intestine. Administration of GH to mice resulted in an increase in GRTP1 mRNA in testes, but a decrease of GRTP1 mRNA in kidney and liver. GRTP1 was identified as an HIV dependency factor (HDF), suggesting that GRTP1 may be an important drug target in HIV treatment. At least three isoforms of GRTP1 are known to exist; this antibody should recognize all isoforms.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	GRTP1 antibody was raised in rabbits against an 18 amino acid peptide from near the amino terminus of human GRTP1.
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, ICC, IF Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. GRTP1 antibody can be used for detection of GRTP1 by Western blot at 1 - 2 µg/mL. Antibody can also be used for immunocytochemistry starting at 2.5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Lu C, Kasik J, Stephan DA, et al. Grtp1, a novel gene regulated by growth hormone. <i>Endocrinology</i> 2001; 142:4568-71. Brass AL, Dykxhoorn DM, Benita Y, et al. Identification of host proteins required for HIV infection through a functional genomic screen. <i>Science</i> 2008; 319:921-6.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4589P; GRTP1 Peptide Cat. No. PK-AB718-1204; K562 Cell Lysate

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