

GPR15 / BOB (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat G-Protein Coupled Receptor 15 (HIV Coreceptor)

Instruction Manual

Catalog Number	PK-AB718-2101
Synonyms	GPR15 Antibody: BOB
Description	Human immunodeficiency virus (HIV) and related viruses require coreceptors to infect target cells. Some G protein-coupled receptors including CCR5, CXCR4, CCR3, CCR2b and CCR8 in the chemokine receptor family, and four new human molecules GPR15, STRL33, GPR1 and V28 were recently identified as HIV coreceptors. The genes encoding human and monkey GPR15/BOB (for G protein-coupled receptor 15 and brother of Bonzo, respectively,) were recently cloned. This novel G protein-coupled receptor serves as coreceptor for simian immunodeficiency virus (SIV), and for strains of HIV-2 and M-tropic HIV-1. The ligand for GPR15 has not been identified yet.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal GPR15 antibody was raised against a peptide corresponding to amino acids 13 to 28 of human GPR15. The sequence differs from those of African green monkey and pig-tailed macaque BOB by one amino acid
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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.h application have to be determined individually. GPR15 antibody can be used for detection of GPR15/BOB by Western blot at 0.5 - 1 µg/mL. An approximately 50 kDa band can be detected.
Images	Available upon request.
References	Heiber M, Marchese A, Nguyen T, Heng HH, George SR, O'Dowd BF. A novel human gene encoding a G-protein-coupled receptor (GPR15) is located on chromosome 3. Genomics 1996;32:462-5 Deng HK, Unutmaz D, KewalRamani VN, Littman DR. Expression cloning of new receptors used by simian and human immunodeficiency viruses. Nature 199;388:296-300 Farzan M, Choe H, Martin K, Marcon L, Hofmann W, Karlsson G, Sun Y, Barrett P, Marchand N, Sullivan N, Gerard N, Gerard C, Sodroski J. Two orphan seven-transmembrane segment receptors which are expressed in CD4-positive cells support simian immunodeficiency virus infection. J Exp Med 1997;186:405-11
Images	Available upon request.
Related Products	CCR5 Antibody, Cat. No. PK-AB718-1112 CCR8 Antibody, Cat. No. PK-AB718-2097 CXCR4 Antibody, Cat. No. PK-AB718-1009 GPR15 Peptide, Cat. No. PK-AB718-2101P Spleen Lysate, Cat. No. PK-AB718-1306

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