

Instruction Manual

Catalog Number	PK-AB718-2209
Synonyms	Bonzo Antibody: STRL33
Description	Human immunodeficiency virus (HIV) and simian immunodeficiency virus (SIV) require coreceptors, in addition to CD4, to infect target cells. Some G protein-coupled receptors including CCR5, CXCR4, CCR3, and CCR2b in the chemokine receptor family have been identified as HIV coreceptors. An orphan G protein-coupled receptor was recently cloned and designated Bonzo, STRL33 and TYMSTR, and identified as HIV and SIV coreceptor. Bonzo/STRL33 serves as coreceptor for SIV, HIV-2 and HIV-1. The messenger RNA of Bonzo/STRL33 is expressed in lymphoid tissues and activated peripheral blood lymphocytes.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Bonzo antibody was raised against a peptide corresponding to amino acids 11 to 25 of human Bonzo/STRL33. The sequence of this peptide differs from those of African green monkey and pig-tailed macaque by one or two amino acids, respectively, .
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.h application have to be determined individually.
Images	Bonzo antibody can be used for detection of Bonzo by Western blot at 0.5 µg/mL. Antibody can also be used for immunohistochemistry starting at 20 µg/mL. For immunofluorescence start at 20 µg/mL.
References	Deng HK, Unutmaz D, KewalRamani VN, Littman DR. Expression cloning of new receptors used by simian and human immunodeficiency viruses. <i>Nature</i> 1997;388:296-300 Liao F, Alkhatib G, Peden KW, Sharma G, Berger EA, Farber JM. STRL33, A novel chemokine receptor-like protein, functions as a fusion cofactor for both macrophage-tropic and T cell line-tropic HIV-1. <i>J Exp Med</i> 1997;185:2015-23 Alkhatib G, Liao F, Berger EA, Farber JM, Peden KW. A new SIV co-receptor, STRL33. <i>Nature</i> 1997;388:238 Loetscher M, Amara A, Oberlin E, Brass N, Legler D, Loetscher P, D'Apuzzo M, Meese E, Rousset D, Virelizier JL, Baggiolini M, Arenzana-Seisdedos F, Moser B. TYMSTR, a putative chemokine receptor selectively expressed in activated T cells, exhibits HIV-1 coreceptor function. <i>Curr Biol</i> 1997;7:652-60
Images	Available upon request.
Related Products	Bonzo Peptide, Cat. No. PK-AB718-2209P; Spleen Lysate, Cat. No. PK-AB718-1306 Bonzo Antibody, Cat. No. PK-AB718-1170; CXCR4 Antibody, Cat. No. PK-AB718-1009 CCR3 Antibody, Cat. No. PK-AB718-1105; CCR5 Antibody, Cat. No. PK-AB718-1112.

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