

# CXCR4-Lo antibody (pAb)

## Rabbit Anti-Human CXCR4-Lo

### Instruction Manual

Catalog Number	PK-AB718-4443
Synonyms	CXCR4-Lo Antibody; CXCR4-Lo, Fusin, LESTR, HUMSTR
Description	Human immunodeficiency virus (HIV) and related viruses require coreceptors, in addition to CD4, 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. Among them, CXCR4 is a principal coreceptor for T-cell tropic strains of HIV-1 fusion and entry of human white blood cells. CXCR4 is also required for the infection by dual-tropic strains of HIV-1 and mediates CD-4 independent infection by HIV-2. The $\alpha$ -chemokine SDF-1 is the ligand for CXCR4 and prevents infection by T-tropic HIV-1. CXCR4 associates with the surface CD4-gp120 complex before HIV enters target cells. CXCR4 messenger RNA levels correlated with HIV-1 permissiveness in diverse human cell types. Antibodies to CXCR4 block HIV-1 and HIV-2 fusion and infection of human target cells. The amino-terminal domain and the second extracellular loop of CXCR4 serve as HIV binding sites. This antibody is specific for the longer isoform of CXCR4.
Quantity	100 $\mu$ g
Source / Host	Rabbit
Immunogen	CXCR4-Lo antibody was raised in rabbits against a peptide corresponding to nine amino acids near the amino terminus of human CXCR4 isoform a.
Purification Method	Purified IgG
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	This antibody is specific for the longer isoform of CXCR4.
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 $\mu$ l 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. CXCR4-Lo antibody can be used for Western blot at 10 $\mu$ g/mL. Antibody can also be used for immunocytochemistry starting at 2 $\mu$ g/mL. For immunofluorescence start at 4 $\mu$ g/mL.
Images	Available upon request.
References	Dimitrov DS. Cell 1997; 91:721-30. Feng Y et al. Science 1996; 272:872-7. Berson JF et al. J. Virol. 1996; 70:6288-95. Doranz BJ et al. Cell 1996; 85:1149-58.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4443P; CXCR4-Lo Peptide Cat.No. PK-AB718-1306; Human Spleen Tissue Lysate

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