

HIV-Igp120 (NT) antibody (pAb)

Rabbit Anti-Glycoprotein 120 (NT)

Instruction Manual

Catalog Number	PK-AB718-4641
Synonyms	gp120 Antibody: HIV-1 glycoprotein 120
Description	Human immunodeficiency virus type 1 (HIV-1) entry into target cells is directed by the envelope (Env) glycoproteins, which are present on the surface of HIV-1 virion or infected cells in the form of trimers consisting of gp120/gp41 complexes. The surface subunit, gp120, initiates the entry process by interacting sequentially with the CD4 receptor and a co-receptor CCR5 or CXCR4, thereby inducing a conformational change that allows the transmembrane (TM) gp41 subunit to mediate fusion between viral and target cell membranes. Cleavage of Env into its gp120 and gp41 components is necessary for activation of its fusogenic activity.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	gp120 antibody was raised in rabbits against 16 amino acid peptide representing the major HIV-1 consensus sequence near the N-terminal region of V1/V2 loop.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Virus
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 Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Gp120 antibody can be used for detection of gp120 by Western blot at 0.5 - 1 µg/mL.
Images	Available upon request.
References	Pinter A. Roles of HIV-1 Env variable regions in viral neutralization and vaccine development. <i>Curr. HIV Res.</i> 2007; 5:542-53. Alkhatib G and Berger EA. HIV coreceptors: from discovery and designation to new paradigms and promise. <i>Eur. J. Med. Res.</i> 2007; 12:375-84.
Images	Available upon request.
Related Products	Cat.No. PK-AB718-4641P; gp120 Peptide

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