

DC-SIGN (CT) antibody (pAb)

Rabbit Anti-Human DC-SIGN (Novel HIV Binding Protein)

Instruction Manual

Catalog Number	PK-AB718-2347
Synonyms	DC-SIGN Antibody:
Description	Dendritic cells (DCs) that control immune responses were recently found to capture and transport HIV from the mucosal area to remote lymph nodes, where DCs hand over HIV to CD4+ T lymphocytes. DCs also amplify the amount of virus and extend the duration of viral infectivity. Multiple strains of HIV-1, HIV-2 and SIV bind to DCs via DC-SIGN. ICAM-3 is the natural ligand for DC-SIGN. A DC-SIGN homologue (termed DC-SIGNR, L-SIGN, and DC-SIGN2) was identified recently. DC-SIGN forms a novel gene family with DC-SIGNR and many alternatively spliced isoforms of DC-SIGN and DC-SIGNR. The expression of DC-SIGN was found in mucosal tissues including placenta, small intestine, and rectum.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal DC-SIGN antibody was raised against a synthetic peptide corresponding to amino acids near the center of human DC-DIGN .
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. DC-SIGN antibody can be used for detection of DC-SIGN by Western blot at 1 to 2 µg/mL. A band at approximately 44 kDa can be detected. Antibody can also be used for immunohistochemistry starting at 10 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Geijtenbeek TB, Kwon DS, Torensma R, et al. DC-SIGN, a dendritic cell-specific HIV-1-binding protein that enhances trans-infection of T cells. <i>Cell</i> . 2000;100:587-97. Pohlmann S, Baribaud F, Lee B, et al. RW. DC-SIGN interactions with human immunodeficiency virus type 1 and 2 and simian immunodeficiency virus. <i>J Virol</i> . 2001;75(10):4664-72. Geijtenbeek TB, Torensma R, van Vliet SJ, et al. Identification of DC-SIGN, a novel dendritic cell-specific ICAM-3 receptor that supports primary immune responses. <i>Cell</i> . 2000;100(5):575-85. Soilleux EJ, Barten R, Trowsdale J. DC-SIGN; a related gene, DC-SIGNR; and CD23 form a cluster on 19p13. <i>J Immunol</i> . 2000;165(6):2937-42.
Images	Available upon request.
Related Products	Blocking Peptide, Cat. No. PK-AB718-2347P Human Small Intesting Tissue Lysate, Cat. No. PK-AB718-1308 DC-SIGN Antibody (ED), Cat. No. PK-AB718-2349

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