

CX3CL1 / CX3C antibody (pAb)

Rabbit Anti-Human/Mouse CX3C-Chemokine Ligand 1 (CX3CR1 Ligand, Neurotactin, Fractalkine)

Instruction Manual

Catalog Number	PK-AB718-2099
Synonyms	CX3CL1 Antibody: CX3C, CX3C membrane-anchored chemokine, neurotactin, small
Description	Chemokines are a family of proteins associated with the trafficking of leukocytes in immune surveillance and inflammatory cell recruitment. They are classified based on the positions of key cysteine residues. CX3CL1 is a CX3C chemokine known to induce adhesion and migration of leukocytes mediated by a membrane-bound and soluble form respectively. Recent experiments have shown that CX3CL1 can suppress the production of nitrous oxide, interleukin-6, and TNF- α in activated microglia and neuronal cells, suggesting that it may act as an intrinsic inhibitor against neurotoxicity by activated microglia. Its receptor, CX3CR1, also functions as a co-receptor for HIV-1 and HIV-2 envelope fusion and virus infection, which can be inhibited by CX3CL1. CX3CL1 is often observed migrating at 80 – 100 kDa in SDS-PAGE, presumably due to post-translational modification.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	CX3CL1 antibody was raised against a peptide corresponding to 18 amino acids near the amino terminus of human CX3CL1.
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse
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 μ 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 INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. CX3CL1 antibody can be used for detection of CX3CL1 by Western blot at 0.5 - 2 mg/mL dilution. CX3CL1 is often observed migrating at 80 - 100 kDa in SDS-PAGE, presumably due to post-translational modification.
Images	Available upon request.
References	Bajetto A, Bonavia R, Barbero S, et al. Chemokines and their receptors in the central nervous system. <i>Front. Neuroendocrinol.</i> 2001; 22:147-84. Umehara H, Goda S, Imai T, et al. Fractalkine, a CX3C-chemokine, functions predominantly as an adhesion molecule in monocytic cell line THP-1. <i>Immunol. Cell. Biol.</i> 2001; 79:298-302. Mizuno T, Kawanokuchi J, Numata K, et al. Production and neuroprotective functions of fractalkine in the central nervous system. <i>Brain Res.</i> 2003; 979:65-70. Combadiere C, Salzwedel K, Smith ED, et al. Identification of CX3CL1. A chemotactic receptor for the human CX3C chemokine fractalkine and a fusion coreceptor for HIV-1. <i>J. Biol. Chem.</i> 1998; 273:23799-804.
Images	Available upon request.
Related Products	C2C12 Lysate, Cat. No. PK-AB718-1285 CX3CL1 Peptide, Cat. No. PK-AB718-2099P

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