

## CX3CR1 Antibody [#G062]

<b>Catalog Number:</b>	G062
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P49238
<b>Other Names:</b>	CX3C chemokine receptor 1; C-X3-C CKR-1; CX3CR1; Fractalkine receptor; G-protein coupled receptor 13; V28; Beta chemokine receptor-like 1; CMK-BRL-1; CMKBLR1; CX3CR1; CMKBRL1; GPR13
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human CX3CR1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CX3CR1 antibody detects endogenous levels of total CX3CR1 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Raport C.J., Gene 163:295-299(1995). Combadiere C., DNA Cell Biol. 14:673-680(1995). DeVries M.E., J. Biol. Chem. 278:11985-11994(2003).