

CCR4 Antibody [#C12128]

Catalog Number:	C12128
Concentration:	1mg/ml
Swiss-Prot No.:	P51679
Other Names:	C-C chemokine receptor type 4, C-C CKR-4, CC-CKR-4, CCR-4, CCR4, K5-5, CD194 antigen, CCR4, CMKBR4
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized peptide derived from internal of human CCR4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CCR4 antibody detects endogenous levels of total CCR4 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 83%, Positives = 100%); Rat (Identities = 83%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:60000
References:	Power C.A., J. Biol. Chem. 270:19495-19500(1995). Kato H., Genes Immun. 1:97-104(1999). Imai T., J. Biol. Chem. 272:15036-15042(1997).