

CCRL2 Antibody [#G061]

Catalog Number:	G061
Concentration:	1mg/ml
Swiss-Prot No.:	O00421
Other Names:	C-C chemokine receptor type 2; C-C CKR-2; CC-CKR-2; CCR-2; CCRL2; CKR2; CMKBR2; JE/FIC receptor; MCP-1 receptor; MCP-1-R; Monocyte chemoattractant protein 1 receptor
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 CCRL2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CCRL2 antibody detects endogenous levels of total CCRL2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:20000
References:	Fan P., Biochem. Biophys. Res. Commun. 243:264-268(1998). Muzny D.M., Nature 440:1194-1198(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).