

## CCR7 Antibody [#G053]

<b>Catalog Number:</b>	G053
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P32248
<b>Other Names:</b>	BLR2; C-C chemokine receptor type 7; C-C CKR-7; CC-CKR-7; CCR-7; CMKBR7; EB11; EBV-induced G-protein coupled receptor 1; MIP-3 beta receptor
<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 CCR7.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	CCR7 antibody detects endogenous levels of total CCR7 protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Birkenbach M.P., J. Virol. 67:2209-2220(1993). Schweickart V.L., Genomics 23:643-650(1994). Imami K., Anal. Sci. 24:161-166(2008).