

IL20RB Antibody [#C12439]

Catalog Number:	C12439
Concentration:	1mg/ml
Swiss-Prot No.:	Q6UXL0
Other Names:	DIRS1; FNDC6; I20RB; IL-20R-beta; IL-20R2; interleukin 20 receptor beta
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 IL20RB.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	IL20RB antibody detects endogenous levels of total IL20RB protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Clark H.F., Genome Res. 13:2265-2270(2003). The MGC Project Team; Genome Res. 14:2121-2127(2004). Zhang Z., Protein Sci. 13:2819-2824(2004).