

## C1QB Antibody [#C15196]

<b>Catalog Number:</b>	C15196
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P02746
<b>Other Names:</b>	complement C1q subcomponent B; complement component 1, q subcomponent, B
<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 C1QB.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	C1QB antibody detects endogenous levels of total C1QB protein.
<b>Reactivity:</b>	Human
<b>Applications:</b>	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Reid K.B.M.; Biochem. J. 231:729-735(1985). Gregory S.G., Nature 441:315-321(2006). The MGC Project Team; Genome Res. 14:2121-2127(2004).