

## ABL1 Antibody [#C10256]

<b>Catalog Number:</b>	C10256
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P00519
<b>Other Names:</b>	ABL1, EC 2.7.1.112, JTK7, Proto-oncogene tyrosine-protein kinase ABL1, c-ABL, p150
<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 ABL1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	ABL1 antibody detects endogenous levels of total ABL1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
<b>References:</b>	Chisoe S.L., Genomics 27:67-82(1995). Wisniewski D., Blood 93:2707-2720(1999). Lin W.-H., Genomics 74:12-20(2001).