

## c-Abl (Ab-412) Antibody [#B7003]

<b>Catalog Number:</b>	B7003
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9NYB9
<b>Other Names:</b>	Abelson interactor 2; ABI2; Abl binding protein 3; Abl-interactor 2; AblBP3; Arg binding protein 1; ArgBP1; ARGBPIA
<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 non-phosphopeptide derived from human c-Abl around the phosphorylation site of tyrosine 412 (D-T-Y <sup>P</sup> -T-A).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	c-Abl (Ab-412) antibody detects endogenous levels of total c-Abl protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
<b>References:</b>	Dai Z., Genes Dev. 9:2569-2582(1995). Ren R.; Submitted (JUL-1995) to the EMBL/GenBank/DDBJ databases. Wang B., Oncogene 12:1921-1929(1996).