

## B-RAF Antibody [#C10378]

<b>Catalog Number:</b>	C10378
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P15056
<b>Other Names:</b>	B-RAF, B-RAF proto-oncogene serine/threonine-protein kinase, B-Raf proto-oncogene serine/threonine-protein kinase, BRAF, BRAF1, EC 2.7.1.37, EC 2.7.11.1, RAFB1, RMIL, RMIL serine/threonine-protein kinase, c-RMIL, kinase B-Raf, p94, v-Raf murine sarcoma viral oncogene homolog B1
<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 B-RAF.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	B-RAF antibody detects endogenous levels of total B-RAF protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Stephens R.M., Mol. Cell. Biol. 12:3733-3742(1992). Hillier L.W., Nature 424:157-164(2003). Eychene A., Oncogene 7:1657-1660(1992).