

B-Raf Antibody [#B0780]

Catalog Number:	B0780
Concentration:	1mg/ml
Swiss-Prot No.:	P15056
Other Names:	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
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 B-Raf.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	B-Raf antibody detects endogenous levels of total B-Raf protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Armelle Calipel, J. Biol. Chem., Oct 2003; 278: 42409 - 42418. Tara J. Dillon, J. Biol. Chem., Sep 2003; 278: 35940 - 35949. R K Jaiswal, Mol. Cell. Biol., Oct 1994; 14: 6944 - 6953. Laura L. Dugan, J. Biol. Chem., Sep 1999; 274: 25842.