

B-RAF (Ab-446) Antibody [#B8082]

Catalog Number:	B8082
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; c-RML; EC 2.7.1.37; EC 2.7.11.1; kinase B-Raf; p94; RAFB1; RML; RML serine/threonine-protein kinase; 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 non-phosphopeptide derived from human B-RAF around the phosphorylation site of serine 446 (R-D-S ^P -S-D).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	B-RAF (Ab-446) antibody detects endogenous levels of total B-RAF protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:40000
References:	Stephens R.M., Mol. Cell. Biol. 12:3733-3742(1992). Albert S., Submitted (JAN-2004) to the EMBL/GenBank/DBJ databases. Hillier L.W., Nature 424:157-164(2003).