

A-RAF (Ab-301/302) Antibody [#B0770]

Catalog Number:	B0770
Concentration:	1mg/ml
Swiss-Prot No.:	P10398
Other Names:	A-RAF, A-Raf proto-oncogene serine/threonine-protein kinase, A-raf-1, ARAF, ARAF1, EC 2.7.11.1, PKS, PKS2, Proto-oncogene Pks, Raf, kinase A-Raf
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 A-RAF around the phosphorylation site of tyrosine 301/tyrosine 302 (G-Y-Y ^P -W-E).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	A-RAF (Ab-301/302) antibody detects endogenous levels of total A-RAF protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
References:	Oliver G. Rössler. Am J Physiol Cell Physiol, May 2004; 286: C1118 - C1129. Marie A. Bogoyevitch. J. Biol. Chem., Nov 1995; 270: 26303. Weicheng Liu. J. Biol. Chem., Mar 2004; 279: 10167 - 10175. M Huleihel. Mol. Cell. Biol., Jul 1986; 6: 2655 - 2662.