

RASSF2 Antibody [#C18180]

Catalog Number:	C18180
Concentration:	1mg/ml
Swiss-Prot No.:	P50749
Other Names:	ras association domain family 2; RASF2; rasfadin
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 RASSF2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RASSF2 antibody detects endogenous levels of total RASSF2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 93%, Positives = 100%); Rat (Identities = 93%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Burbee D.G., Submitted (SEP-2002) to the EMBL/GenBank/DDBJ databases. Nagase T., DNA Res. 3:17-24(1996). Ota T., Nat. Genet. 36:40-45(2004).