

RASSF4 Antibody [#C18182]

Catalog Number:	C18182
Concentration:	1mg/ml
Swiss-Prot No.:	Q9H2L5
Other Names:	RAB5 interacting protein 2; Ras and Rab interactor 2; RAS association (RalGDS/AF-6) domain containing protein JC265; Ras association domain family 4; Ras inhibitor JC265; Ras interaction/interference protein 2; RASSF4
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 RASSF4.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RASSF4 antibody detects endogenous levels of total RASSF4 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Burbee D.G., Submitted (JAN-2003) to the EMBL/GenBank/DDBJ databases. Xu X., Submitted (MAY-2000) to the EMBL/GenBank/DDBJ databases. Deloukas P., Nature 429:375-381(2004).