

RHBT2 Antibody [#C11647]

Catalog Number:	C11647
Concentration:	1mg/ml
Swiss-Prot No.:	Q9BYZ6
Other Names:	Rho-related BTB domain-containing protein 2, Deleted in breast cancer 2 gene protein p83
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 RHBT2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	RHBT2 antibody detects endogenous levels of total RHBT2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:5000
References:	Ramos S., Gene 298:147-157(2002). Nagase T., DNA Res. 5:277-286(1998). Nakajima D., DNA Res. 9:99-106(2002).