

GRB2 Antibody [#C15972]

Catalog Number:	C15972
Concentration:	1mg/ml
Swiss-Prot No.:	P62993
Other Names:	Growth factor receptor-bound protein 2; Adapter protein GRB2; SH2/SH3 adapter GRB2; Protein Ash
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 GRB2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GRB2 antibody detects endogenous levels of total GRB2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Lowenstein E.J., Cell 70:431-442(1992). Matuoka K., Proc. Natl. Acad. Sci. U.S.A. 89:9015-9019(1992). Fath I., Science 264:971-974(1994).