

GRAP2 Antibody [#C15966]

Catalog Number:	C15966
Concentration:	1mg/ml
Swiss-Prot No.:	O75791
Other Names:	GRB2-related adapter protein 2, GADS protein; Growth factor receptor-binding protein; GRBLG; Grf40 adapter protein; Grf-40; GRB-2-like protein; GRB2L; GRBX; P38; Hematopoietic cell-associated adapter protein GrpL; Adapter protein GRID; SH3-SH2-SH3 adapter Mona
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 GRAP2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GRAP2 antibody detects endogenous levels of total GRAP2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Qiu M., Biochem. Biophys. Res. Commun. 253:443-447(1998). Law C.-L., J. Exp. Med. 189:1243-1253(1999). Asada H., J. Exp. Med. 189:1383-1390(1999).