

G3BP2 Antibody [#C18193]

Catalog Number:	C18193
Concentration:	1mg/ml
Swiss-Prot No.:	Q9UN86
Other Names:	G3B2; G3BP2; GAP SH3-domain binding protein 2; KIAA0660; Ras-GTPase-activating protein binding protein 2
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 C-terminal of human G3BP2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	G3BP2 antibody detects endogenous levels of total G3BP2 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Kennedy D., Submitted (APR-1999) to the EMBL/GenBank/DDBJ databases. Ishikawa K., DNA Res. 5:169-176(1998). Kennedy D., Submitted (MAR-1998) to the EMBL/GenBank/DDBJ databases.