

GPRIN1 Antibody [#C15835]

Catalog Number:	C15835
Concentration:	1mg/ml
Swiss-Prot No.:	Q7Z2K8
Other Names:	G protein-regulated inducer of neurite outgrowth 1; KIAA1893
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 GPRIN1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	GPRIN1 antibody detects endogenous levels of total GPRIN1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Nagase T., DNA Res. 8:179-187(2001). The MGC Project Team; Genome Res. 14:2121-2127(2004). The German cDNA consortium; Submitted (JUL-2002) to the EMBL/GenBank/DDBJ databases.