

CG028 Antibody [#G223]

Catalog Number:	G223
Concentration:	1mg/ml
Swiss-Prot No.:	O95766
Other Names:	UPF0550 protein C7orf28; C7orf28A; CGI-43; C7orf28B
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 CG028.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CG028 antibody detects endogenous levels of total CG028 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Lai C.-H., Genome Res. 10:703-713(2000). Ota T., Nat. Genet. 36:40-45(2004). Hillier L.W., Nature 424:157-164(2003).