

FZD2 Antibody [#G107]

Catalog Number:	G107
Concentration:	1mg/ml
Swiss-Prot No.:	Q14332
Other Names:	Frizzled-2 [Precursor]; Fz-2; hFz2; FzE2; FZD2
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 FZD2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FZD2 antibody detects endogenous levels of total FZD2 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
References:	Zhao Z., Genomics 27:370-373(1995). Sagara N., Biochem. Biophys. Res. Commun. 252:117-122(1998). The MGC Project Team; Genome Res. 14:2121-2127(2004).