

LRP11 Antibody [#C16508]

Catalog Number:	C16508
Concentration:	1mg/ml
Swiss-Prot No.:	Q86VZ4
Other Names:	low density lipoprotein receptor-related protein 11; MANSC3
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 LRP11.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LRP11 antibody detects endogenous levels of total LRP11 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Ota T., Nat. Genet. 36:40-45(2004). Mungall A.J., Nature 425:805-811(2003). The MGC Project Team; Genome Res. 14:2121-2127(2004).