

LRP10 Antibody [#C16507]

Catalog Number:	C16507
Concentration:	1mg/ml
Swiss-Prot No.:	Q7Z4F1
Other Names:	Low-density lipoprotein receptor-related protein 10
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 LRP10.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LRP10 antibody detects endogenous levels of total LRP10 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:10000
References:	Clark H.F., Genome Res. 13:2265-2270(2003). Li W.B., Submitted (JAN-2003) to the EMBL/GenBank/DDBJ databases. Bechtel S., BMC Genomics 8:399-399(2007).