

CEP97 Antibody [#C16453]

Catalog Number:	C16453
Concentration:	1mg/ml
Swiss-Prot No.:	Q8IW35
Other Names:	LRRIQ2; leucine-rich repeat and IQ domain-containing 2
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 CEP97.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	CEP97 antibody detects endogenous levels of total CEP97 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 85%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	The MGC Project Team, Genome Res. 14:2121-2127(2004). Ota T., Nat. Genet. 36:40-45(2004). Spektor A., Cell 130:678-690(2007).