

KLHL29 Antibody [#C16417]

Catalog Number:	C16417
Concentration:	1mg/ml
Swiss-Prot No.:	Q96CT2
Other Names:	Kelch-like protein 29, Kelch repeat and BTB domain-containing protein 9
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 KLHL29.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	KLHL29 antibody detects endogenous levels of total KLHL29 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 92%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Bechtel S., BMC Genomics 8:399-399(2007). The MGC Project Team; Genome Res. 14:2121-2127(2004). Nagase T., DNA Res. 8:179-187(2001).