

LAMA5 Antibody [#C13068]

Catalog Number:	C13068
Concentration:	1mg/ml
Swiss-Prot No.:	O15230
Other Names:	Laminin subunit alpha-5; KIAA0533; KIAA1907
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 LAMA5.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LAMA5 antibody detects endogenous levels of total LAMA5 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:40000
References:	Deloukas P., ure 414:865-871(2001). Nagase T., DNA Res. 8:179-187(2001). Nagase T., DNA Res. 5:31-39(1998).