

LAMA1 Antibody [#C13064]

Catalog Number:	C13064
Concentration:	1mg/ml
Swiss-Prot No.:	P25391
Other Names:	LAMA, laminin A, laminin alpha-1
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 LAMA1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LAMA1 antibody detects endogenous levels of total LAMA1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:40000
References:	Haaparanta T., Matrix 11:151-160(1991). Nissinen M., Biochem. J. 276:369-379(1991). Olsen D., Lab. Invest. 60:772-782(1989).