

## LAMA2 Antibody [#C13065]

<b>Catalog Number:</b>	C13065
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P24043
<b>Other Names:</b>	LAMM, laminin M, laminin alpha-2, merosin heavy chain
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from internal of human LAMA2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	LAMA2 antibody detects endogenous levels of total LAMA2 protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 85%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
<b>References:</b>	Vuolteenaho R., J. Cell Biol. 124:381-394(1994). Ehrig K., . Natl. Acad. Sci. U.S.A. 87:3264-3268(1990). Liu T., J. Proteome Res. 4:2070-2080(2005).