

## LAMC1 Antibody [#C13072]

<b>Catalog Number:</b>	C13072
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P11047
<b>Other Names:</b>	Laminin subunit gamma-1 [Precursor], Laminin B2 chain, LAMC1, LAMB2
<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 LAMC1.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	LAMC1 antibody detects endogenous levels of total LAMC1 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Kallunki T., J. Biol. Chem. 266:221-228(1991). Pikkarainen T., J. Biol. Chem. 263:6751-6758(1988). Fukushima Y., Cytogenet. Cell Genet. 48:137-141(1988).