

LAMB3 Antibody [#C13071]

Catalog Number:	C13071
Concentration:	1mg/ml
Swiss-Prot No.:	Q13751
Other Names:	Laminin subunit beta-3, Laminin 5 beta 3, Laminin B1k chain, Kalinin B1 chain, LAMB3, LAMNB1
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 LAMB3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	LAMB3 antibody detects endogenous levels of total LAMB3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
References:	Gerecke D.R., J. Biol. Chem. 269:11073-11080(1994). Morishima Y., J. Biol. Chem. 270:273-279(1995). Robbins P.B., Proc. Natl. Acad. Sci. U.S.A. 98:5193-5198(2001).