

Collagen IV α 2 Antibody [#C12196]

Catalog Number:	C12196
Concentration:	1mg/ml
Swiss-Prot No.:	P08572
Other Names:	CO4A2, Collagen alpha-2(IV) chain, alpha 2 type IV collagen, canstatin, collagen type IV alpha 2
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 Collagen IV α 2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen IV α 2 antibody detects endogenous levels of total Collagen IV α 2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Dunham A., Nature 428:522-528(2004). Fischer G., Biochem. J. 292:687-695(1993). Kamphaus G.D., J. Biol. Chem. 275:1209-1215(2000).