

Collagen XI α 1 Antibody [#C12214]

Catalog Number:	C12214
Concentration:	1mg/ml
Swiss-Prot No.:	P12107
Other Names:	Collagen alpha-1(XI) chain [Precursor], COBA1
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 XI α 1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen XI α 1 antibody detects endogenous levels of total Collagen XI α 1 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 ELISA: 1:10000
References:	Yoshioka H., J. Biol. Chem. 265:6423-6426(1990). Annunen S., Am. J. Hum. Genet. 65:974-983(1999). Zhidkova N.I., J. Biol. Chem. 270:9486-9493(1995).