

Collagen XIII α 1 Antibody [#C12217]

Catalog Number:	C12217
Concentration:	1mg/ml
Swiss-Prot No.:	Q5TAT6
Other Names:	Collagen alpha-1(XIII) chain, COLXIII A1, CODA1
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 XIII α 1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen XIII α 1 antibody detects endogenous levels of total Collagen XIII α 1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IF: 1:500~1:1000 ELISA: 1:20000
References:	Snellman A., EMBO J. 19:5051-5059(2000). Deloukas P., Nature 429:375-381(2004). Pihlajaniemi T., J. Biol. Chem. 265:16922-16928(1990).