

Collagen I α 2 Antibody [#C12195]

Catalog Number:	C12195
Concentration:	1mg/ml
Swiss-Prot No.:	P08123
Other Names:	Collagen alpha-2(I) chain [Precursor], Alpha-2 type I collagen, CO1A2
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 I α 2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen I α 2 antibody detects endogenous levels of total Collagen I α 2 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:20000
References:	Dalgleish R., Nucleic Acids Res. 25:181-187(1997). Korkko J.M., Am. J. Hum. Genet. 62:98-110(1998). Akai J., Connect. Tissue Res. 30:1-6(1998).