

## Collagen I Antibody [#C0154]

<b>Catalog Number:</b>	C0154
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P02452
<b>Other Names:</b>	CO1A1, alpha 1 type I collagen preproprotein, collagen alpha 1(I) chain, collagen type I alpha 1
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized peptide derived from N-terminal of human Collagen I.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	Collagen I antibody detects endogenous levels of total Collagen I protein.
<b>Reactivity:</b>	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
<b>Applications:</b>	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	Masatoshi Jinnin, Nucleic Acids Res., Mar 2005; 33: 1337 - 1351. B. Stefanovic, Nucleic Acids Res., Jan 2000; 28: 641 - 647. DE Birk, J. Cell Biol., Mar 1988; 106: 999. Christina Meyer, J. Biol. Chem., Nov 2002; 277: 45400 - 45407.