

Collagen XVIII α 1 Antibody [#C12223]

Catalog Number:	C12223
Concentration:	1mg/ml
Swiss-Prot No.:	P39060
Other Names:	Collagen alpha-1(XVIII) chain, Endostatin, COIA1
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 XVIII α 1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen XVIII α 1 antibody detects endogenous levels of total Collagen XVIII α 1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 92%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Saarela J., Matrix Biol. 16:319-328(1998). Elamaa H., Matrix Biol. 22:427-442(2003). Hattori M, Nature 405:311-319(2000).