

Collagen VI α 2 Antibody [#C12205]

Catalog Number:	C12205
Concentration:	1mg/ml
Swiss-Prot No.:	P12110
Other Names:	alpha 2 type VI collagen, collagen alpha-2
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 VI α 2.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen VI α 2 antibody detects endogenous levels of total Collagen VI α 2 protein.
Reactivity:	Human, Mouse
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
References:	The MGC Project Team, Genome Res. 14:2121-2127(2004). Saitta B., J. Biol. Chem. 267:6188-6196(1992). Tillet E., Eur. J. Biochem. 221:177-185(1994).