

Collagen VI α 1 Antibody [#C12204]

Catalog Number:	C12204
Concentration:	1mg/ml
Swiss-Prot No.:	P12109
Other Names:	Collagen alpha-1(VI) chain, CO6A1
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 1.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen VI α 1 antibody detects endogenous levels of total Collagen VI α 1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 81%, Positives = 100%)
Applications:	IHC: 1:50~1:100 ELISA: 1:20000
References:	Tillet E., Eur. J. Biochem. 221:177-185(1994). Trikkka D., Mamm. Genome 8:342-345(1997). Scacheri P.C., Neurology 58:593-602(2002).