

Collagen VI α 3 Antibody [#C12206]

Catalog Number:	C12206
Concentration:	1mg/ml
Swiss-Prot No.:	P12111
Other Names:	Collagen alpha-3(VI) chain, CO6A3
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 3.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Collagen VI α 3 antibody detects endogenous levels of total Collagen VI α 3 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%)
Applications:	IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Chu M.-L., EMBO J. 9:385-393(1990). Zanussi S., J. Biol. Chem. 267:24082-24089(1992). Liu T., J. Proteome Res. 4:2070-2080(2005).