

## Collagen IV Antibody [#C0157]

- Catalog Number:** C0157
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P02462
- All Names:** COL4A1 NC1 domain, COLLAGEN OF BASEMENT MEMBRANE, ALPHA-1 CHAIN ARRESTEN, alpha 1 type IV collagen preproprotein, collagen IV alpha-1 polypeptide, collagen type IV alpha 1
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized peptide derived from human Collagen IV.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Collagen IV antibody detects endogenous levels of total Collagen IV protein.
- Reactivity:** Human, Mouse
- Applications:** WB: 1:300~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
- References:** Mary K. Chelberg, Cancer Res., Sep 1989; 49: 4796 - 4802.  
Patricia L. Graham, J. Cell Biol., Jun 1997; 137: 1171.  
PJ Lein, J. Cell Biol., Apr 1991; 113: 417.  
EC Tsilibary, J. Biol. Chem., Dec 1988; 263: 19112 - 19118.