

## TIMP2 Antibody [#C0348]

- Catalog Number:** C0348-1, C0348-2, C0348
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P16035
- All Names:** tissue inhibitor of metalloproteinase 2
- 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 TIMP2.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** TIMP2 antibody detects endogenous levels of total TIMP2 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IF: 1:500~1:1000 ELISA: 1:40000
- References:** Dominique Trudel, AACR Meeting Abstracts, Mar 2004; 2004: 264.  
O. Alcazar, Invest. Ophthalmol. Vis. Sci., May 2007; 48: 5057.  
Isabelle Kazes, Blood, Nov 2000; 96: 3064 - 3069.  
Georgina S. Butler, J. Biol. Chem., Jan 1998; 273: 871.