

TGF β 1 Antibody [#C0340]

- Catalog Number:** C0340-1, C0340-2, C0340
Amount: 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
Swiss-Prot No. : P01137
All Names: CED, DPD1, TGF-beta 1, TGFB, diaphyseal dysplasia 1, progressive, transforming growth factor, beta 1, transforming growth factor beta 1 (Camurati-Engelmann disease)
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), 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 TGF β 1 antibody.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** TGF β 1 antibody detects endogenous levels of total TGF β 1 protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
- References:** Shu Wu, Am. J. Respir. Cell Mol. Biol., May 2007; 36: 552 - 561.
Chun-Lin Chen, J. Biol. Chem., Apr 2006; 281: 11506 - 11514.
Ying Dong Xu, Am J Physiol Lung Cell Mol Physiol, Sep 2003; 285: 527.
Q H Song, Mol. Pathol., Jun 2002; 55: 164 - 176.