

TGF β 2 Antibody [#C0341]

- Catalog Number:** C0341-1, C0341-2, C0341
Amount: 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
Swiss-Prot No. : P61812
All Names: TRANSFORMING GROWTH FACTOR BETA-2
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 2.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity TGF β 2 antibody detects endogenous levels of total TGF β 2 protein.
Reactivity: Human (Identities = 100%, Positives = 100%);
Mouse (Identities = 95%, Positives = 95%);
Rat (Identities = 95%, Positives = 95%)
Applications: IHC: 1:50~1:100 ELISA: 1:2000
References: C Gordon-Thomson, Invest. Ophthalmol. Vis. Sci., Jul 1998; 39: 1399.
Elisabetta del Re, J. Biol. Chem., May 2004; 279: 22765 - 22772.
H. Garrett R. Thompson, Am J Physiol Lung Cell Mol Physiol, Dec 2006; 291: L1277 - L1285.
FRANCIS J. KELLY, J. Am. Soc. Nephrol., Jun 1999; 10: 1264.