

TGF β Receptor II Antibody [#C0338]

- Catalog Number:** C0338-1, C0338-2, C0338
Amount: 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
Swiss-Prot No. : P37173
All Names: GF-beta type II receptor, TGF-betaR2, TGFR-2, TGFR2, kinase TGF-beta-R2
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 Receptor II antibody.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: TGF β Receptor II antibody detects endogenous levels of total TGF β Receptor II protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:20000
References: Herbert Y. Lin, J. Biol. Chem., Feb 1995; 270: 2747.
Michitomo Sakuma, Int. Immunol., Feb 2007; 19: 117 - 126.
Elisabetta del Re, J. Biol. Chem., May 2004; 279: 22765 - 22772.
M Laiho, J. Biol. Chem., Oct 1990; 265: 18518 - 18524.