

FGF Receptor 1 (Ab-154) Antibody [#B7084]

- Catalog Number:** B7084-1, B7084-2, B7084
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : P11362
All Names: EC 2.7.10.1, FGFBR, FGFR-1, FGR1, FLG, FLT2, Fms-like tyrosine kinase-2, MFR, bFGF-R, c-fgr, kinase FGFR1
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 non-phosphopeptide derived from human FGF Receptor 1 around the phosphorylation site of tyrosine 154 (A-P-Y-W-T).
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity FGF Receptor 1 (Ab-154) antibody detects endogenous levels of total FGF Receptor 1 protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:40000
References: Vikash Kumar Dubey, Protein Sci., Sep 2005; 14: 2315 - 2323.
F. Mouriaux, Invest. Ophthalmol. Vis. Sci., May 2004; 45: 1206.
E Shi, Mol. Cell. Biol., Jul 1993; 13: 3907 - 3918.
Dara J. Dunican, Cell Growth Differ., May 2001; 12: 255.