

FGFR1 Oncogene Partner Antibody [#C0190]

Catalog Number: C0190-1, C0190-2, C0190

Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl

Swiss-Prot No.: O95684

All Names: FOP, FIBROBLAST GROWTH FACTOR RECEPTOR 1 ONCOGENE PARTNER

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 FGFR1 Oncogene Partner.

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity FGFR1 Oncogene Partner antibody detects endogenous levels of total FGFR1 Oncogene Partner protein.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000

References: Asuman Demiroglu, Blood, Dec 2001; 98: 3778 - 3783.

Géraldine Guasch, Blood, Jan 2003; 101: 286 - 288.

Ronald Simon, Cancer Res., Jun 2001; 61: 4514 - 4519.

Ute Schell, Hum. Mol. Genet., Mar 1995; 4: 323 - 328.