

PDGF Receptor β (Ab-752) Antibody [#B7194]

Catalog Number:	B7194
Concentration:	1mg/ml
Swiss-Prot No.:	P09619
Other Names:	CD140b; EC 2.7.10.1; kinase PDGFR-beta; PDGF-R-beta; PDGFR; PDGFR-beta; PGFRB; platelet-derived growth factor receptor, beta polypeptide
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg^{2+} and Ca^{2+}), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human PDGF Receptor β around the phosphorylation site of tyrosine 751 (V-D-Y ^P -V-P).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PDGF Receptor β (Ab-752) antibody detects endogenous levels of total PDGF Receptor β protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Gronwald R.G.K., Proc. Natl. Acad. Sci. U.S.A. 85:3435-3439(1988). Claesson-Welsh L., Mol. Cell. Biol. 8:3476-3486(1988). The MGC Project Team; Genome Res. 14:2121-2127(2004).