

## PDGFR $\beta$ (Ab-740) Antibody [#B0970]

<b>Catalog Number:</b>	B0970
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P09619
<b>Other Names:</b>	CD140b, EC 2.7.10.1, PDGF-R-beta, PDGFR, PDGFR-beta, PGFRB, kinase PDGFR-beta
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human PDGFR $\beta$ around the phosphorylation site of tyrosine 740 (G-G-Y <sup>P</sup> -M-D).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	PDGFR $\beta$ (Ab-740) antibody detects endogenous levels of total PDGFR $\beta$ protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	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). Chi K.D., Oncogene 15:1051-1057(1997).