

PDGFRa (Ab-849) Antibody [#B8210]

Catalog Number:	B8210
Concentration:	1mg/ml
Swiss-Prot No.:	P16234
Other Names:	Alpha platelet-derived growth factor receptor precursor; CD140a; EC 2.7.10.1; kinase PDGFR-alpha; PDGF-R-alpha; PDGFR-alpha; PGFRA; platelet-derived growth factor, alpha-receptor
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 PDGFRa around the phosphorylation site of tyrosine 849 (S-N-Y ^P -V-S).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PDGFRa (Ab-849) antibody detects endogenous levels of total PDGFRa protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Mouse (Identities = 100%, Positives = 100%); Rat (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Claesson-Welsh L., Proc. Natl. Acad. Sci. U.S.A. 86:4917-4921(1989). Matsui T., Science 243:800-804(1989). Kawagishi J., Genomics 30:224-232(1995).