

PDGFRa(Phospho-Tyr849)Antibody [#8A8210]

Catalog Number:	8A8210
Concentration:	1mg/ml
Origin of Antibody:	Rabbit
Swiss-Prot No:	P16234
NCBI Gene Symbol:	PGFRA
Other Names:	Alpha platelet-derived growth factor receptor precursor; CD140a; kinase PDGFR-alpha; PDGF-R-alpha; PDGFR-alpha; PGFRA; platelet-derived growth factor, alpha-receptor
Modification Sites:	H:Y849 M:Y849 R:Y848
Target Modification:	Phospho
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 peptide derived from human PDGFRa around the phosphorylation site of Tyr849.
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
Specificity:	PDGFRa (Phospho-Tyr849) Antibody detects endogenous levels of PDGFRa only when phosphorylated at Tyr849.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 ELISA: 1:1000