

## Data Sheet

### PDGFR $\alpha$

Human, recombinant, N-terminal GST tag

**Catalog #:** 40261

Lot #: 110811

Conc.: 0.1mg/ml

**Formulated in:** 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.1 mM EGTA, 0.1 mM PMSF, and 25% glycerol, and 0.25 mM DTT.

**Stability:** >6 months at  $-80^{\circ}\text{C}$

**References:**

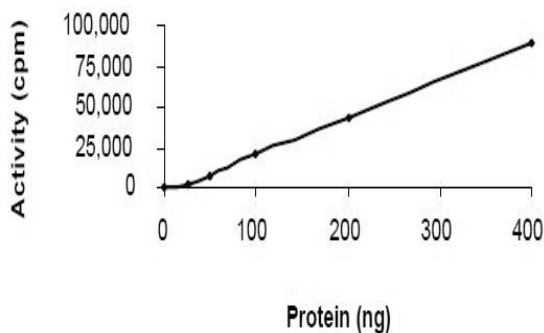
1. Floege, J. *et al. J Am Soc Nephrol.* 1998 Feb; **9(2)**:211-23.
2. Kumagai, S. *et al. Tohoku J Exp Med.* 2001 Sep; **195(1)**:21-33.

**Description:** Recombinant human PDGFR $\alpha$  (550-end), Genbank accession number NM\_00626, with an N-terminal GST tag, expressed by baculovirus in Sf9 insect cells.

**Specific Activity:** 12 pmol/min/ $\mu\text{g}$ .  
 Assay condition: 25 mM MOPS, 12.5 mM  $\beta$ -glycerophosphate, 20 mM  $\text{MgCl}_2$ , 25 mM  $\text{MnCl}_2$ , 5 mM EGTA, 2 mM EDTA, 0.25 mM DTT, 50  $\mu\text{g/ml}$  BSA, 2 mM ATP, PDGFR $\alpha$ , and 1 mg/ml poly Glu:Tyr (4:1) peptide as substrate. Incubate with [ $^{33}\text{P}$ ]-ATP at  $30^{\circ}\text{C}$  for 15 minutes, then spot on phosphocellulose paper, fix in 1% phosphoric acid, and assay with a scintillation counter.

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance



**10% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
 PDGFR $\alpha$

**Lane 2:**  
 Protein Marker

**MW:** 95 kDa  
**Purity:**  $\geq 90\%$

