

Data Sheet

PDGFR α

Mouse, recombinant, N-terminal GST tag

Catalog #: 40260

Lot #: 110519

Conc.: 0.1mg/ml

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

Stability: >6 months at -80°C

References:

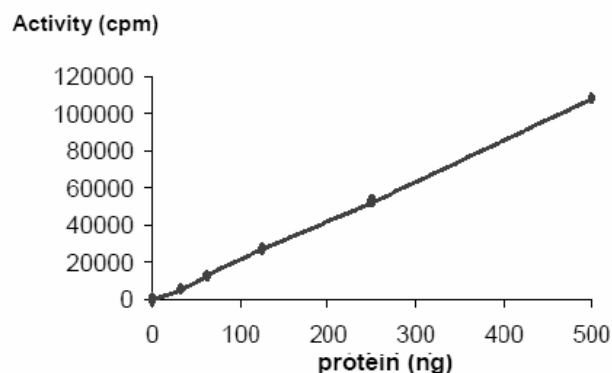
1. Floege, J. *et al. J Am Soc Nephrol.* 1998 Feb; **9(2)**:211-23.
2. Kumagai, S. *et al. Tohoku J Exp Med.* 2001 Sep; **195(1)**:21-33.

Description: Recombinant murine PDGFR α (550-end), Genbank accession number NM_011058, with an N-terminal GST tag, expressed by baculovirus in Sf9 insect cells.

Specific Activity: 14 pmol/min/ μg .
 Assay condition: 5 mM MOPS, 2.5 mM β -glycerophosphate, 5 mM MgCl_2 , 1 mM EGTA, 0.4 mM EDTA, 0.05 mM DTT, 50 $\mu\text{g/ml}$ BSA, 2 mM ATP, PDGFR α , and 1 mg/ml poly Glu:Tyr (4:1) peptide as substrate. Incubate with [^{33}P]-ATP at 30°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 α 5 μg

Lane 2:
 Protein Marker

MW: 100 kDa
Purity: $\geq 90\%$

