

## Data Sheet

### PDGFR $\beta$

Human, recombinant, N-terminal His-tag

**Catalog #:** 40263

**Lot#:** 100812 **Conc.:** 0.66 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM Imidazole, 3 mM DTT, and 10% glycerol.

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

**References:**

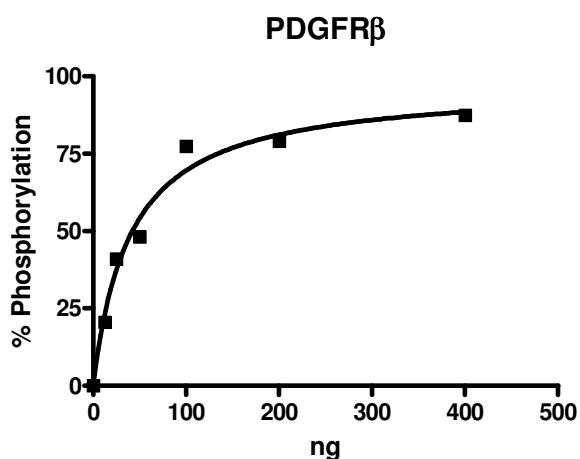
1. Claesson-Welsh L. *et al.* (1988). *Mol. Cell. Biol.* **8**:3476.
2. Heldin C.-H. and Westermark B. (1999). *Physiol. Rev.* **79**:1283.

**Description:** Human PDGFR $\beta$  (GenBank Accession No. NM\_002609), cytoplasmic domain, a.a. 558-1106(end) with N-terminal His-tag, MW= 62 kDa, expressed in Sf9 cells via a Baculovirus expression system.

**Specific Activity:** 15 pmol/min/ $\mu\text{g}$ .  
**Assay condition:** The enzyme reaction was carried out for 1h at  $30^{\circ}\text{C}$  in a buffer containing 50 mM HEPES (pH7.5), 10 mM  $\text{MgCl}_2$ , 1 mM EDTA, 0.01% BRIJ-35 and 200  $\mu\text{M}$  of ATP. Substrate peptide: Tyr2 peptide from Invitrogen.

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

### **Quality Assurance**



**10% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
 PDGFR $\beta$  5  $\mu\text{g}$

**Lane 2:**  
 Protein Marker  
 BioLabs (#P7708L)

**MW:** 62 kDa.  
**Purity:**  $\geq 50\%$

