

PDGFR β

(PGFRB, Beta platelet-derived growth factor receptor [Precursor])

- **CAT#:** PDGFR β
- **Enzyme:** Human PDGFR β
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	85
PP2	30,000
PDGFR I-III	3,300
D-64406	1,600

- **Scientific Information:** P09619 (<http://www.expasy.org/uniprot/P09619>)

Reference Compound IC50 for PDGFR β

