

PDGFR α

(PGFRA, Alpha platelet-derived growth factor receptor)

- **CAT#:** PDGFR α
- **Enzyme:** Human PDGFR α
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 10 μ M
- **Reaction:**

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

Compounds	IC50 (nM)
Staurosporine	< 0.1
PP2	18.9
AG1478	306
PDGFR I-III	2.4
D-64406	453

- **Scientific Information:** Q9DE49 (<http://www.expasy.org/uniprot/Q9DE49>)

Reference compound IC50 for PDGFR α

