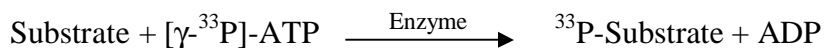


PDGFR α

(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 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.2
PP2	900
AG1478	9,700
PDGFR I-III	614
D-64406	310

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

Reference Compound IC50 for PDGFR α

