

PDGFR α (V561D)

(PGFRA, Alpha platelet-derived growth factor receptor)

- **CAT#:** PDGFR α (V561D)
- **Enzyme:** Human PDGFR α (V561D)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 10 μ M
- **Reaction:**

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

Compounds	IC50 (nM)
Staurosporine	< 0.1
PP2	12
AG1478	328
PDGFR I-III	1.9
D-64406	217

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

Reference compound IC50 for PDGFR α (V561D)

