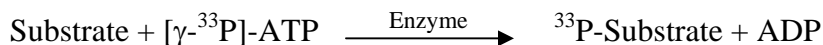


## PDGFRa (T674I)

(PGFRA, Alpha platelet-derived growth factor receptor)

- **CAT#:** PDGFRa (T674I)
- **Enzyme:** Human PDGFRa (T674I)
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.8
PP2	>50,000
AG1478	821
PDGFR I-III	67
D-64406	2,600

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

### Reference compound IC50 for PDGFRa (T674I)

