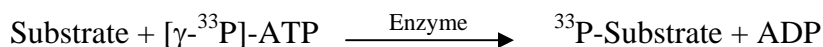


## TEC

(Tyrosine-protein kinase TEC, PSCTK4)

- **CAT#:** TEC
- **Enzyme:** Human TEC
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	600
PP2	4,500
D-64406	18,000

- **Scientific Information:** P42680 (<http://www.expasy.org/uniprot/P42680>)

### Reference compound IC50 for TEC

