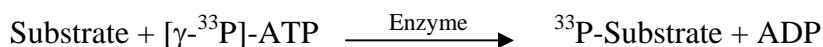


## TXK

(Tyrosine-protein kinase TXK, PTK4)

- **CAT#:** TXK
- **Enzyme:** Human TXK
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	11.7
PP2	3.6
AG1478	923
D-64406	4,300

- **Scientific Information:** P42681 (<http://www.expasy.org/uniprot/P42681>)

### Reference compound IC50 for TXK

