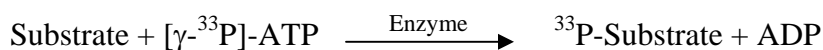


## TYRO3/SKY

(Tyrosine-protein kinase receptor TYRO3 [Precursor], Tyrosine-protein kinase RSE, Tyrosine-protein kinase SKY, Tyrosine-protein kinase DTK, Protein-tyrosine kinase byk, BYK, DTK, RSE, SKY)

- **CAT#:** TYRO3
- **Enzyme:** Human TYRO3
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.1
PP2	5,800
AG1478	2,900
D-64406	3,700

- **Scientific Information:** Q06418 (<http://www.expasy.org/uniprot/Q06418>)

**Reference Compound IC50 for TYRO3/SKY**

