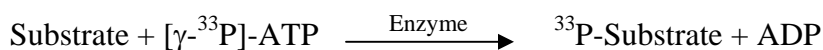


TYRO3/SKY

(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/SKY
- **Enzyme:** Human TYRO3/SKY
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	5.6
PP2	>20,000
AG1478	5,100
D-64406	3,200

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

Reference compound IC50 for TYRO3/SKY

