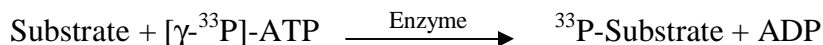


## EPHA2

(Ephrin type-A receptor 2 [Precursor], Tyrosine-protein kinase receptor ECK, Epithelial cell kinase)

- **CAT#:** EPHA2
- **Enzyme:** Human EPHA2
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	274
PP2	99
AG1478	2,400
D-64406	27,000

- **Scientific Information:** P29317 (<http://www.expasy.org/uniprot/P29317>)

### Reference Compound IC50 for EPHA2

