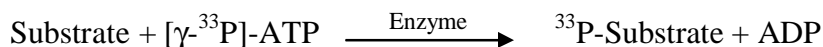


EPHA1

(Ephrin type-A receptor 1 [Precursor], Tyrosine-protein kinase receptor EPH)

- **CAT#:** EPHA1
- **Enzyme:** Human EPHA1
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	74
PP2	21.5
AG1478	2,800
D-64406	34,000

- **Scientific Information:** P21709 (<http://www.expasy.org/uniprot/P21709>)

Reference Compound IC50 for EPHA1

