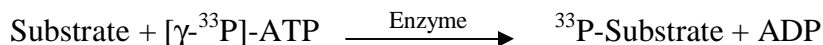


EphB1

(Ephrin type-B receptor 1 [Precursor], Tyrosine-protein kinase receptor EPH-2, NET, HEK6, ELK)

- **CAT#:** EphB1
- **Enzyme:** Human EphB1
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	190
PP2	70
AG1478	756
D-64406	8,800

- **Scientific Information:** P54762 (<http://www.expasy.org/uniprot/P54762>)

Reference Compound IC50 for EPHB1

