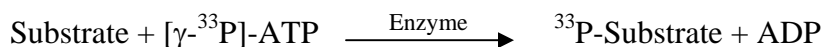


EPHB4

(Ephrin type-B receptor 4 [Precursor], Tyrosine-protein kinase receptor HTK, TYRO11)

- **CAT#:** EPHB4
- **Enzyme:** Human EPHB4
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	112
PP2	96
AG1478	700
D-64406	13,000

- **Scientific Information:** P54760 (<http://www.expasy.org/uniprot/P54760>)

Reference compound IC50 for EPHB4

