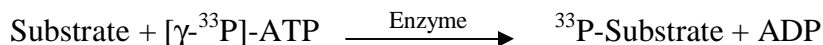


EphB2

(Ephrin type-B receptor 2 [Precursor], Tyrosine-protein kinase receptor EPH-3, DRT, Receptor protein-tyrosine kinase HEK5, ERK, NY-REN-47 antigen)

- **CAT#:** EphB2
- **Enzyme:** Human EphB2
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	30
PP2	13.5
AG1478	473
D-64406	2,600

- **Scientific Information:** P29323 (<http://www.expasy.org/uniprot/P29323>)

Reference Compound IC50 for EPHB2

