

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), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	92
PP2	169
AG1478	2,600
D-64406	22,700

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

Reference compound IC50 for EPHB2

