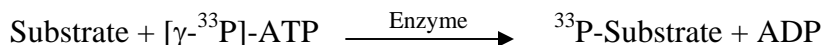


ErbB4/HER4

(Receptor tyrosine-protein kinase erbB-4 [Precursor], p180erbB4,
Tyrosine kinase-type cell surface receptor HER4)

- **CAT#:** ErbB4
- **Enzyme:** Human ErbB4
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	30.9
PP2	62.1
AG1478	26
D-64406	5,900

- **Scientific Information:** Q15303 (<http://www.expasy.org/uniprot/Q15303>)

Reference compound IC50 for ErbB4/HER4

