

## 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 10  $\mu$ M
- **Reaction:**  

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	96.7
PP2	274
AG1478	29.6
D-64406	15,000

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

### Reference compound IC50 for ErbB4/HER4

