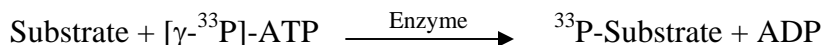


HER2/ERBB2

(Receptor tyrosine-protein kinase erbB-2 [Precursor], p185erbB2, C-erbB-2, NEU proto-oncogene, Tyrosine kinase-type cell surface receptor HER2, MLN 19)

- **CAT#:** HER2/ERBB2
- **Enzyme:** Human HER2/ERBB2
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	106
PP2	214
AG1478	14.5

- **Scientific Information:** P04626 (<http://www.expasy.org/uniprot/P04626>)

Reference Compound IC50 for HER2

