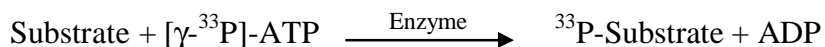


EGFR

(Epidermal growth factor receptor, Receptor tyrosine-protein kinase ErbB-1)

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

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 38 |
| PP2 | 29 |
| AG1478 | 4.5 |
| D-64406 | 6,900 |

- **Scientific Information:** P00533 (<http://www.expasy.org/uniprot/P00533>)

Reference Compound IC50 for EGFR

