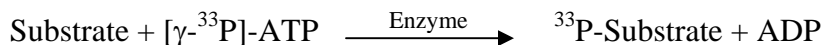


EGFR (L858R, T790M)

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

- **CAT#:** EGFR (L858R, T790M)
- **Enzyme:** Human EGFR (L858R, T790M)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.6
PP2	3,800
AG1478	4,500

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

Reference compound IC50 for EGFR (L858R, T790M)

