

## EGFR (d746-750)

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

- **CAT#:** EGFR (d746-750)
- **Enzyme:** Human EGFR (d746-750)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	27
PP2	320
AG1478	0.2
PDGFR I-III	>50,000
D-64406	6,300

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

### Reference compound IC50 for EGFR (d746-750)

