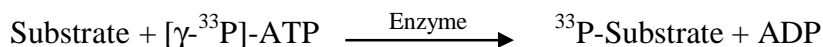


FGFR2

(Fibroblast growth factor receptor 2 [Precursor], Keratinocyte growth factor receptor 2)

- **CAT#:** FGFR2
- **Enzyme:** Human FGFR2
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.4
PP2	940
AG1478	2,500
PDGFR I-III	16,400
D-64406	1,400

- **Scientific Information:** P21802 (<http://www.expasy.org/uniprot/P21802>)

Reference Compound IC50 for FGFR2

