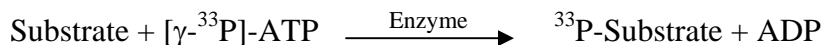


FGFR4

(Fibroblast growth factor receptor 4 [Precursor], FGFR-4, JTK2, TKF)

- **CAT#:** FGFR4
- **Enzyme:** Human FGFR4
- **Substrate:** Peptide substrate, , poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	110
PP2	2,400
AG1478	>20,000
D-64406	>20,000

- **Scientific Information:** P22455 (<http://www.expasy.org/uniprot/P22455>)

Reference compound IC50 for FGFR4

