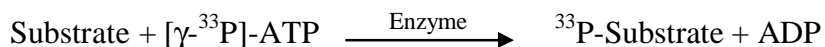


FGFR3

(Fibroblast growth factor receptor 3 [Precursor], FGFR-3)

- **CAT#:** FGFR3
- **Enzyme:** Human FGFR3
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	52
PP2	10,800
D-64406	5,900

- **Scientific Information:** P22607 (<http://www.expasy.org/uniprot/P22607>)

Reference Compound IC50 for FGFR3

