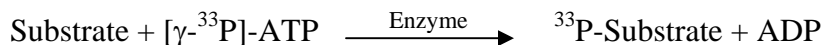


## FGFR3 (K650E)

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

- **CAT#:** FGFR3 (K650E)
- **Enzyme:** Human FGFR3 (K650E)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 5.1       |
| PP2           | 6,100     |
| AG1478        | 3,700     |
| D-64406       | 11,200    |

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

### Reference compound IC50 for FGFR3 (K650E)

