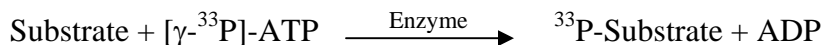


FGFR2 (N549H)

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

- **CAT#:** FGFR2 (N549H)
- **Enzyme:** Human FGFR2 (N549H)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 1 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.4
PP2	61.7
AG1478	307
PDGFR I-III	16,700
D-64406	729

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

Reference Compound IC50 for FGFR2 (N549H)

