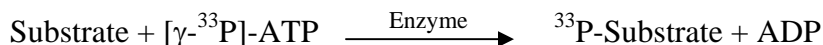


FGFR1 (V561M)

(Basic fibroblast growth factor receptor 1 [Precursor], FGFR-1, bFGF-R, Fms-like tyrosine kinase 2, FLT2, c-fgr)

- **CAT#:** FGFR1 (V561M)
- **Enzyme:** Human FGFR1 (V561M)
- **Substrate:** Peptide substrate, [GEEEEYFELVKKKK], 20 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.1
PDGFR I-III	9,400
D-64406	1,400

- **Scientific Information:** P11362 (<http://www.expasy.org/uniprot/P11362>)

Reference Compound IC50 for FGFR1 (V561M)

