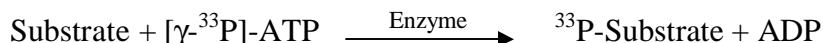


FGFR1

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

- **CAT#:** FGFR1
- **Enzyme:** Human FGFR1
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.3
PP2	1,200
AG1478	12,000
PDGFR I-III	275,000
D-64406	5,700

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

Reference Compound IC50 for FGFR1

