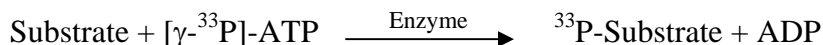


FLT3

(FL cytokine receptor [Precursor], Tyrosine-protein kinase receptor FLT3, Stem cell tyrosine kinase 1, STK-1, Flk-2)

- **CAT#:** FLT3
- **Enzyme:** Human FLT3
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.2
PP2	1,500
AG1478	709
PDGFR I-III	278
D-64406	48.4

- **Scientific Information:** P36888 (<http://www.expasy.org/uniprot/P36888>)

Reference compound IC50 for FLT3

