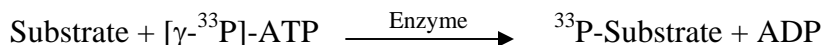


FLT4/VEGFR3

(Vascular endothelial growth factor receptor 3 [Precursor], VEGFR-3,
Tyrosine-protein kinase receptor FLT4, VGFR3)

- **CAT#:** FLT4
- **Enzyme:** Human FLT4
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.2
PP2	900
AG1478	2,200
PDGFR I-III	15,600
D-64406	2,600

- **Scientific Information:** P35916 (<http://www.expasy.org/uniprot/P35916>)

Reference compound IC50 for FLT4/VEGFR3

