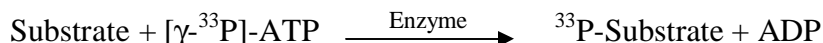


FLT1/VEGFR-1

(Vascular endothelial growth factor receptor 1 [Precursor], Vascular permeability factor receptor, Tyrosine-protein kinase receptor FLT, Tyrosine-protein kinase FRT, Flt-1, Fms-like tyrosine kinase 1)

- **CAT#:** FLT1
- **Enzyme:** Human FLT1
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.5
PP2	1,500
AG1478	2,400
PDGFR I-III	2,700
D-64406	700

- **Scientific Information:** P17948 (<http://www.expasy.org/uniprot/P17948>)

Reference Compound IC50 for FLT1

