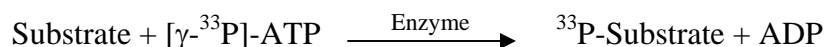


KDR/VEGFR-2

(Vascular endothelial growth factor receptor 2 [Precursor], Kinase insert domain receptor, Protein-tyrosine kinase receptor Flk-1, VEGFR2)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.3
PP2	83.8
AG1478	84.9
PDGFR I-III	4,000
D-64406	454

- **Scientific Information:** P35968 (<http://www.expasy.org/uniprot/P35968>)

Reference compound IC50 for KDR/VEGFR2

