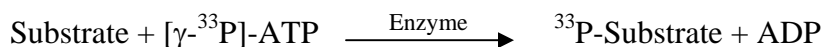


c-Kit (V654A)

(Mast/stem cell growth factor receptor, Proto-oncogene tyrosine-protein kinase c-kit)

- **CAT#:** c-Kit (V654A)
- **Enzyme:** Human c-Kit (V654)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	12
PP2	3,800
AG1478	14,600
PDGFR I-III	2,100
D-64406	144

- **Scientific Information:** P10721 (<http://www.expasy.org/uniprot/P10721>)

Reference Compound IC50 for c-Kit (V654A)

