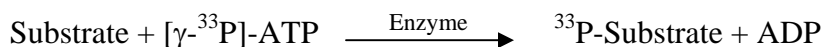


c-Kit (V560G)

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

- **CAT#:** c-Kit (V560G)
- **Enzyme:** Human c-Kit (V560G)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 1 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.14
PP2	50
AG1478	2,000
PDGFR I-III	4.4
D-64406	180

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

Reference Compound IC50 for c-KIT (V560G)

