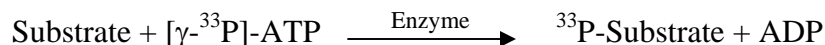


c-Kit

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

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4.2
PP2	114
AG1478	1,800
PDGFR I-III	31
D-64406	61.3

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

Reference compound IC50 for c-KIT

