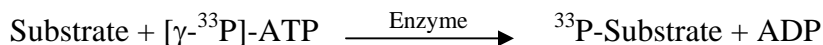


RET

(Proto-oncogene tyrosine-protein kinase receptor ret [Precursor], C-ret)

- **CAT#:** RET
- **Enzyme:** Human RET
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.2
PP2	21
AG1478	350
PDGFR I-III	>20,000
D-64406	250

- **Scientific Information:** P07949 (<http://www.expasy.org/uniprot/P07949>)

Reference compound IC50 for RET

