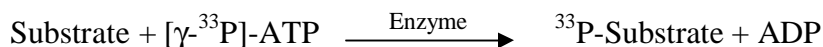


## RET

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

- **CAT#:** RET
- **Enzyme:** Human RET
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 5  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	9.9
PP2	34.8
AG1478	580
D-64406	441

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

**Reference Compound IC50 for RET**

