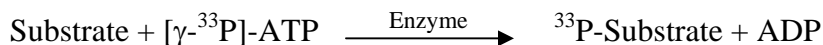


RET (V804M)

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

- **CAT#:** RET (V804M)
- **Enzyme:** Human RET (V804M)
- **Substrate:** Peptide substrate, [KKKVSRSGLYRSPS MPENLNRPR], 20 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.5
PP2	5,000
AG1478	23,700
PDGFR I-III	13,500
D-64406	260

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

Reference Compound IC50 for RET (V804M)

