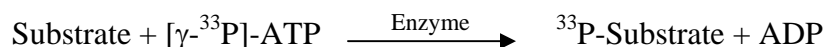


c-MET

(Hepatocyte growth factor receptor [Precursor], HGF receptor, Scatter factor receptor, SF receptor, HGF/SF receptor, Met proto-oncogene tyrosine kinase, c-Met, MET)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	260
PP2	>20,000
AG1478	4,800
PDGFR I-III	>20,000
D-64406	13,000

- **Scientific Information:** P08581 (<http://www.expasy.org/uniprot/P08581>)

Reference compound IC50 for c-MET

