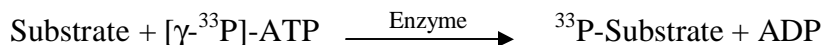


## c-MET

(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], 5  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	216
PP2	52,500
AG1478	51,000
PDGFR I-III	46,000
D-64406	20,600

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

### Reference Compound IC50 for c-MET

