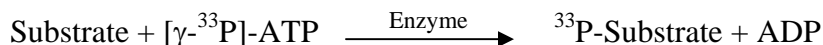


c-MET (M1250T)

(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 (M1250T)
- **Enzyme:** Human c-MET (M1250T)
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	338
AG1478	45,500
D-64406	26,500

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

Reference compound IC50 for c-MET (M1250T)

