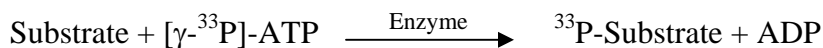


## c-MER

(Proto-oncogene tyrosine-protein kinase MER [Precursor], Receptor tyrosine kinase MerTK, MERTK, MER)

- **CAT#:** c-MER
- **Enzyme:** Human c-MER
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4.5
PP2	3,200
AG1478	2,600
PDGFR I-III	807
D-64406	6,800

- **Scientific Information:** Q12866 (<http://www.expasy.org/uniprot/Q12866>)

**Reference Compound IC50 for c-MER**

