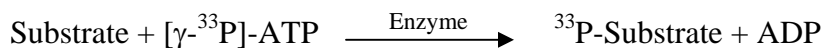


c-MER

(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 10 μ M

- **Reaction:**

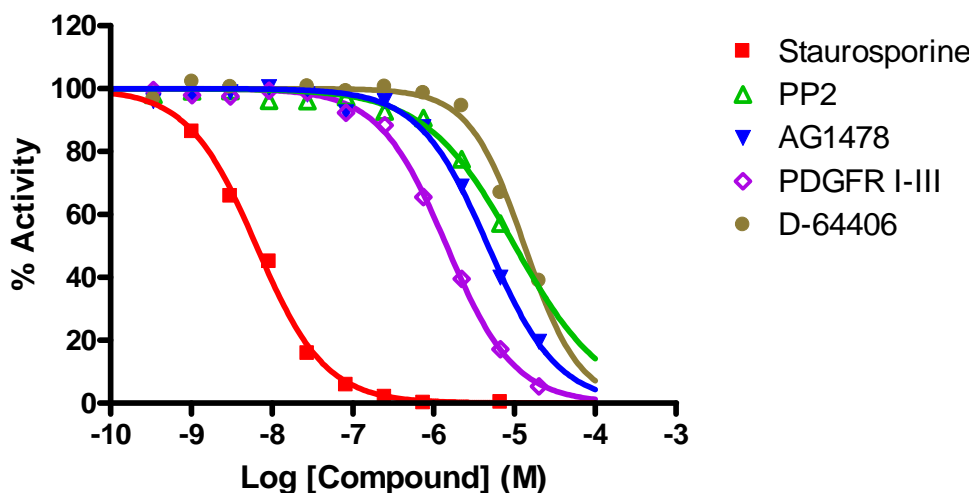


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6.4
PP2	10,100
AG1478	4,700
PDGFR I-III	1,400
D-64406	13,300

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

Reference compound IC50 for c-MER



Service performed
using kinases
under license from  **invitrogen**[™]
part of *life* technologies[™]