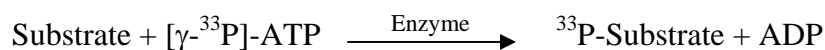


MEK1

(Dual specificity mitogen-activated protein kinase kinase 1, MAP kinase kinase 1, MAPKK 1, ERK activator kinase 1, MAPK/ERK kinase 1, MP2K1, MAP2K1, PRKMK1)

- **CAT#:** MEK1
- **Enzyme:** Human MEK1
- **Substrate:** ERK2 (K52R), 1 μ M
ATP 10 μ M

- **Reaction:**

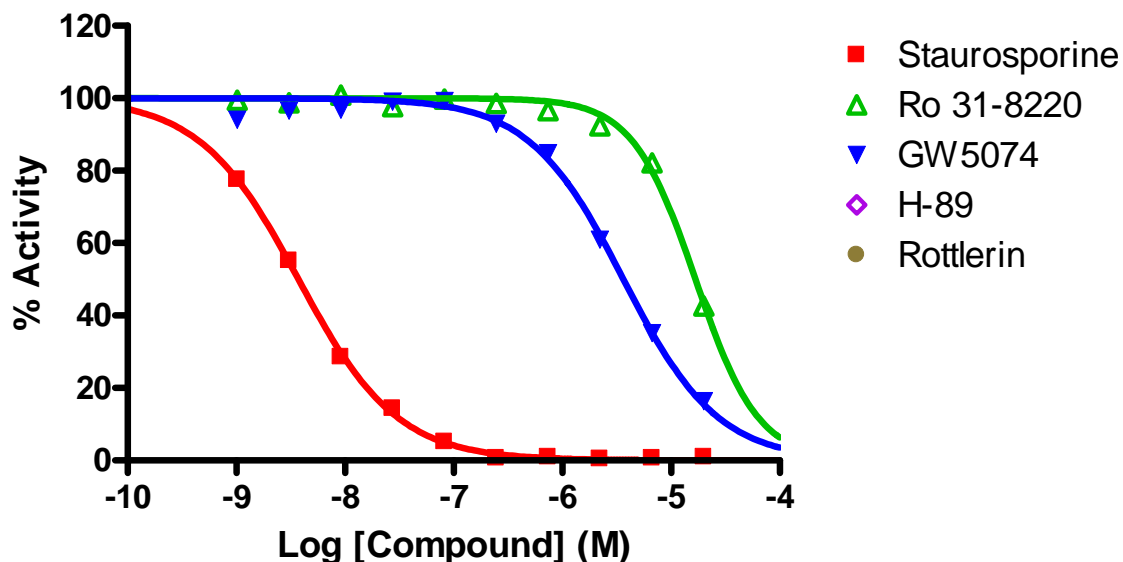


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.7
Ro 31-8220	16,800
GW 5074	3,600

- **Scientific Information:** Q02750 (<http://www.expasy.org/uniprot/Q02750>)

Reference compound IC50 for MEK1



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