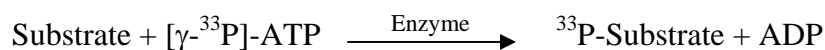


MEK1

(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 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.52
Ro 31-8220	8,700
GW 5074	1,300
H-89	16,700

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

Reference Compound IC50 for MEK1

