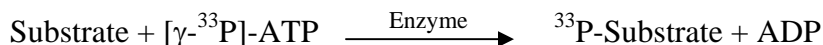


MEKK1

(Mitogen-activated protein kinase kinase kinase 1, MAPK/ERK kinase kinase 1, MEK kinase 1, MAP3K1, MAPKKK1, M3K1)

- **CAT#:** MEKK1
- **Enzyme:** Human MEKK1
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1,200
Ro 31-8220	>20,000
GW 5074	>20,000
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** Q13233 (<http://www.expasy.org/uniprot/Q13233>)

Reference compound IC50 for MEKK1

