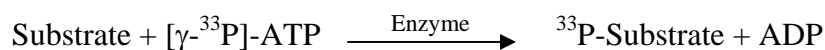


ASK1/MAP3K5

(Mitogen-activated protein kinase kinase kinase 5, MAPK/ERK kinase kinase 5, MEK kinase 5, MEKK 5, Apoptosis signal-regulating kinase 1, ASK-1, M3K5, MAP3K5, MAPKKK5)

- **CAT#:** ASK1/MAP3K5
- **Enzyme:** Human ASK1/MAP3K5
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	24
Ro 31-8220	21,000
GW 5074	23,000
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** Q99683 (<http://www.expasy.org/uniprot/Q99683>)

Reference compound IC50 for ASK1/MAP3K5

