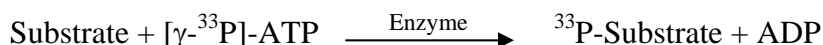


KHS/MAP4K5

(Mitogen-activated protein kinase kinase kinase kinase 5, MAPK/ERK kinase kinase kinase 5, MEK kinase kinase 5, MEKKK 5, Kinase homologous to SPS1/STE20, MAP4K5, M4K5)

- **CAT#:** KHS/MAP4K5
- **Enzyme:** Human KHS/MAP4K5
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.6
Ro 31-8220	260
GW 5074	7,600
H-89	38,000
Rottlerin	68,000

- **Scientific Information:** Q9Y4K4 (<http://www.expasy.org/uniprot/Q9Y4K4>)

Reference Compound IC50 for KHS/MAP4K5

