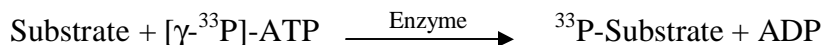


MAPKAPK2

(MAP kinase-activated protein kinase 2, MAPK-activated protein kinase 2, MAPKAP kinase 2, MAPKAPK-2, MK2, MAPK2)

- **CAT#:** MAPKAPK2
- **Enzyme:** Human MAPKAPK2
- **Substrate:** Peptide substrate, [KKLNRTLVA], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	14.5
Ro 31-8220	6,600
GW5074	3,700
H-89	80,000
Rottlerin	9,200

- **Scientific Information:** P49137 (<http://www.expasy.org/uniprot/P49137>)

Reference Compound IC50 for MAPKAPK2

