

MAPKAPK5/PRAK

(MAP kinase-activated protein kinase 5, MAPK-activated protein kinase 5, MAPKAP kinase 5, p38-regulated/activated protein kinase, MAPK5)

- **CAT#:** MAPKAPK5/PRAK
- **Enzyme:** Human MAPKAPK5/PRAK
- **Substrate:** Peptide substrate, [KKLNRTLVA], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	420
Ro 31-8220	15,000
GW 5074	13,000
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** Q8IW41 (<http://www.expasy.org/uniprot/Q8IW41>)

Reference compound IC50 for MAPKAPK5/PRAK

