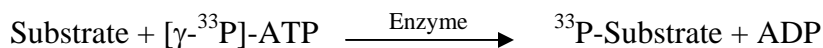


## GCK

(Mitogen-activated protein kinase kinase kinase 2, MAPK/ERK kinase kinase 2, MEK kinase kinase 2, MEKKK 2, Germinal center kinase, GC kinase, Rab8-interacting protein, B lymphocyte serine/threonine-protein kinase, MAP4K2, RAB8IP, M4K2)

- **CAT#:** GCK
- **Enzyme:** Human GCK
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.9
Ro 31-8220	886
GW 5074	2,800
H-89	22,500

- **Scientific Information:** Q12851 (<http://www.expasy.org/uniprot/Q12851>)

Reference compound IC50 for GCK

