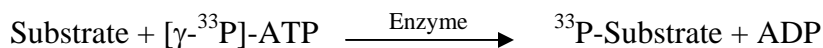


## HGK/MAP4K4

(Mitogen-activated protein kinase kinase kinase kinase 4, MAPK/ERK kinase kinase kinase 4, MEK kinase kinase 4, MEKKK 4, HPK/GCK-like kinase HGK Nck-interacting kinase, M4K4)

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

- **Reaction:**

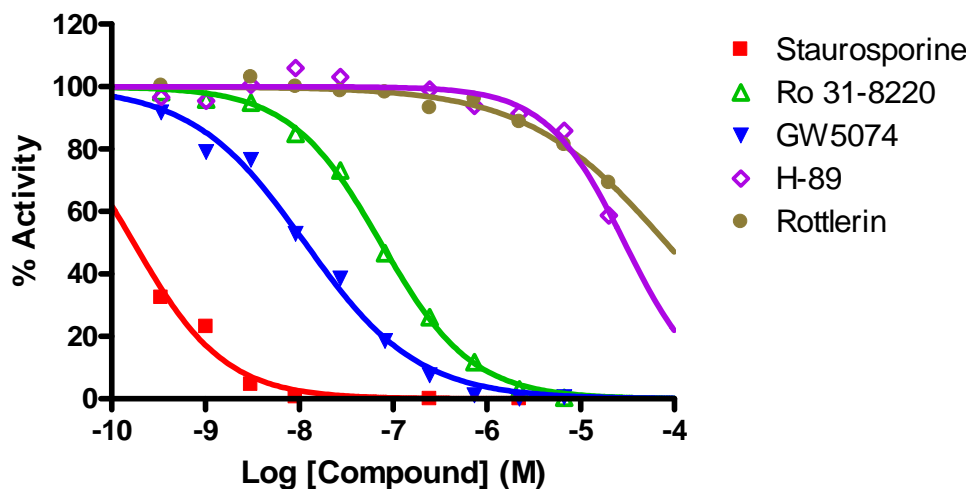


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.17
Ro 31-8220	74.3
GW 5074	11.5
H-89	29,500
Rottlerin	8,200

- **Scientific Information:** O95819 (<http://www.expasy.org/uniprot/O95819>)

### Reference compound IC50s for HGK/MAP4K4



Service performed  
using kinases  
under license from  **invitrogen**<sup>™</sup>  
part of *life* technologies<sup>™</sup>