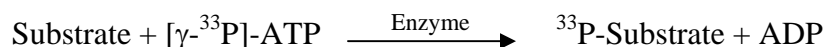


## DLK/MAP3K12

(Mitogen-activated protein kinase kinase kinase 12, Dual leucine zipper bearing kinase, Leucine-zipper protein kinase, MAPK-upstream kinase, Mixed lineage kinase, ZPK, M3K12)

- **CAT#:** DLK/MAP3K12
- **Enzyme:** Human DLK/MAP3K12
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	21
Ro 31-8220	24,000
GW 5074	21,000
H-89	35,000
Rottlerin	>100,000

- **Scientific Information:** Q12852 (<http://www.expasy.org/uniprot/Q12852>)

### Reference compound IC50 for DLK/MAP3K12

