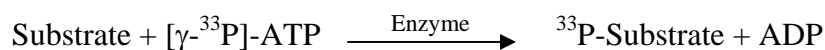


## MEK2

(Dual specificity mitogen-activated protein kinase kinase 2, MAP kinase kinase 2, MAPKK 2, ERK activator kinase 2, MAPK/ERK kinase 2, MP2K2, MAP2K2, PRKMK2, MKK2)

- **CAT#:** MEK2
- **Enzyme:** Human MEK2
- **Substrate:** ERK2 (K52R), 1  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	9.6
Ro 31-8220	8,500
GW 5074	3,900

- **Scientific Information:** P36507 (<http://www.expasy.org/uniprot/P36507>)

### Reference compound IC50 for MEK2

