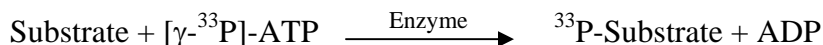


## WNK3

(Serine/threonine-protein kinase WNK3, Protein kinase with no lysine 3  
Protein kinase, lysine-deficient 3)

- **CAT#:** WNK3
- **Enzyme:** Human WNK3
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4,400
Ro 31-8220	>20,000
Wee1 Inhibitor	2,600

- **Scientific Information:** Q9BYP7 (<http://www.expasy.org/uniprot/Q9BYP7>)

### Reference compound IC50 for WNK3

