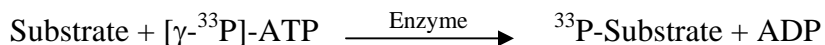


## LCK2/ICK

(Serine/threonine-protein kinase ICK, MAK-related kinase, KIAA0936)

- **CAT#:** LCK2/ICK
- **Enzyme:** Human LCK2/ICK
- **Substrate:** MBP (Myelin Basic Protein), 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	61
Ro 31-8220	40
GW 5074	27,000
H-89	>100,000
Rottlerin	>100,000

- **Scientific Information:** Q9UPZ9 (<http://www.expasy.org/uniprot/Q9UPZ9>)

### Reference compound IC50 for LCK2/ICK

