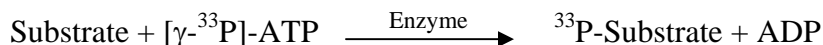


## ZIPK/DAPK3

(Death-associated protein kinase 3, DAP kinase 3, DAP-like kinase  
Dlk, ZIP-kinase)

- **CAT#:** ZIPK
- **Enzyme:** Human ZIPK
- **Substrate:** Peptide substrate, [KKLNRTL<sup>S</sup>FAEPG], 20 μM  
ATP 10 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	5.3
Ro 31-8220	4,000
GW 5074	830
H-89	20,000
Rottlerin	46,000

- **Scientific Information:** O43293 (<http://www.expasy.org/uniprot/O43293>)

### Reference compound IC50 for ZIPK/DAPK3

