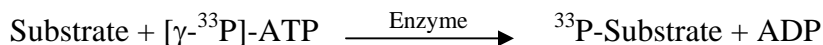


## PKA

(cAMP-dependent protein kinase, PRKACA)

- **CAT#:** PKA
- **Enzyme:** Human PKA
- **Substrate:** Peptide substrate, [GRTGRRNSI-NH<sub>2</sub>], 20 μM  
ATP 10 μM

- **Reaction:**

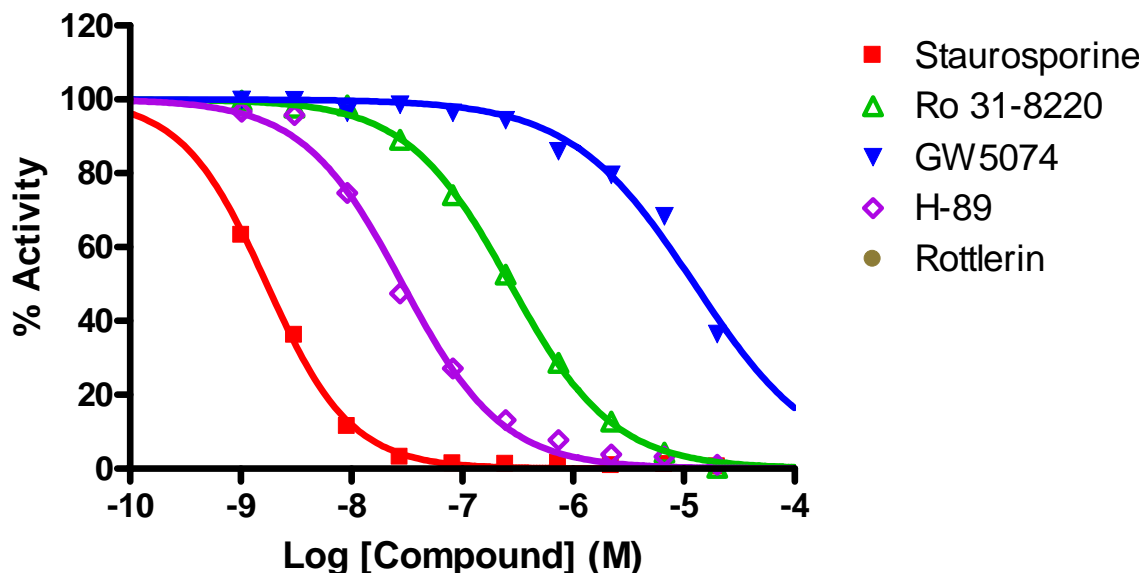


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.7
Ro 31-8220	270
GW 5074	12,000
H-89	29

- **Scientific Information:** P17612 (<http://www.expasy.org/uniprot/P17612>)

### Reference compound IC50 for PKA



Service performed  
using kinases  
under license from  invitrogen™  
part of *Life* technologies™