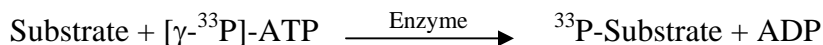


PKC γ

(Protein kinase C gamma type, PKC-gamma, KPCG, PRKCG)

- **CAT#:** PKC γ
- **Enzyme:** Human PKC γ
- **Substrate:** Histone H1, Lipid activator, 20 μ M
ATP 10 μ M

- **Reaction:**

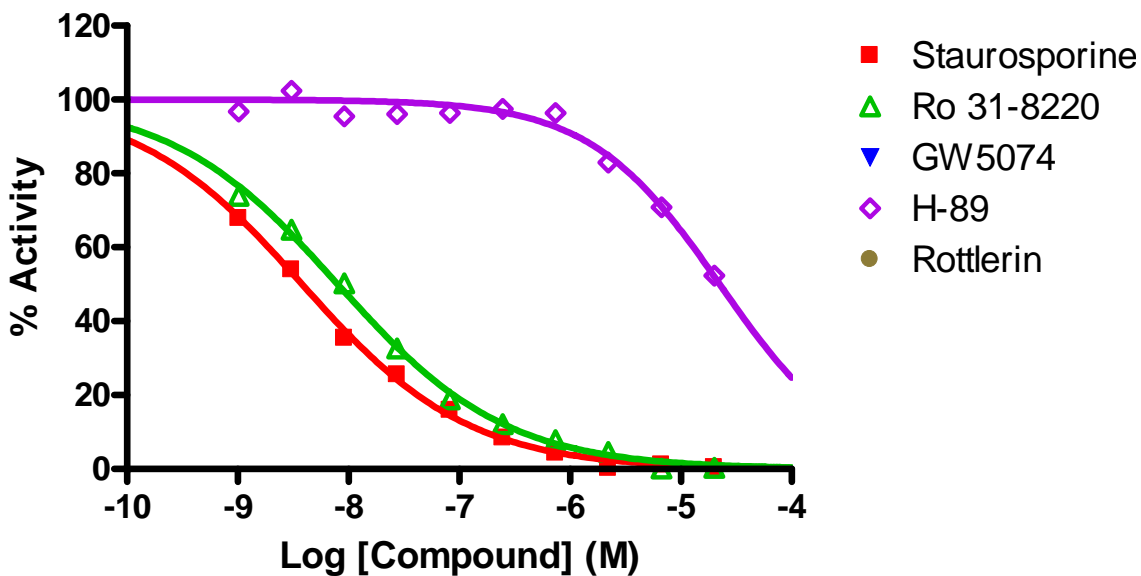


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.8
Ro 31-8220	8.0
H-89	22,000

- **Scientific Information:** P05129 (<http://www.expasy.org/uniprot/P05129>)

Reference compound IC50 for PKC γ



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