

PKC α

(Protein kinase C alpha type, PKC-alpha, PKC-A, KPCA)

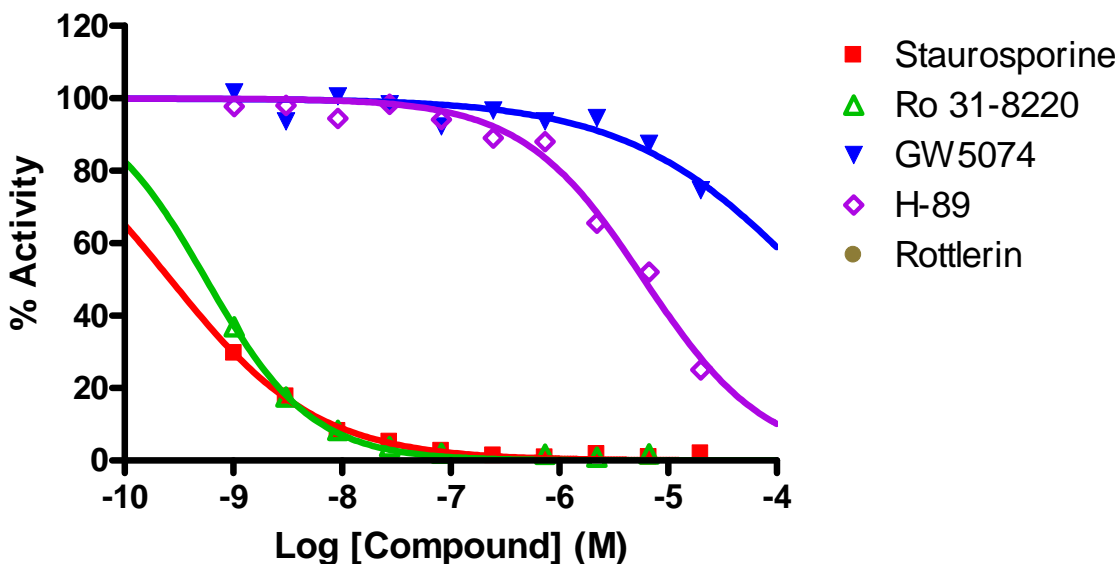
- **CAT#:** PKC α
- **Enzyme:** Human PKC α
- **Substrate:** Histone H1, Lipid activator, 20 μ M
ATP 10 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.26
Ro 31-8220	0.26
GW 5074	>100,000
H-89	6,000

- **Scientific Information:** P17252 (<http://www.expasy.org/uniprot/P17252>)

Reference compound IC50 for PKC α



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