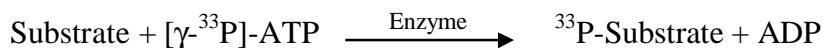


PKC α

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

- **CAT#:** PKC α
- **Enzyme:** Human PKC α
- **Substrate:** Histone H1, Lipid activator, 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.38
Ro 31-8220	0.44
GW 5074	12,000
H-89	2,300
H-89	2,300

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

Reference Compound IC50 for PKC α

