

PKC γ

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

- **CAT#:** PKC γ
- **Enzyme:** Human PKC γ
- **Substrate:** Histone H1, Lipid activator, 5 μ M
ATP 1 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	0.35
Ro 31-8220	0.81
GW5074	7,400
H-89	19,200

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

