

PKC β I

(Protein kinase C beta type, PKC-beta, PKC-B, PRKCB1, KPCB)

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

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

Compounds	IC50 (nM)
Staurosporine	0.75
Ro 31-8220	2.3

- **Scientific Information:** P05771 (<http://www.expasy.org/uniprot/P05771>)

