

## PKC $\eta$

(Protein kinase C eta type, nPKC-eta, PKC-L, KPCL, PRKCH)

- **CAT#:** PKC $\eta$
- **Enzyme:** Human PKC $\eta$
- **Substrate:** Histone H1, Lipid activator, 5  $\mu$ M  
ATP 1  $\mu$ M
- **Reaction:**  

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

Compounds	IC50 (nM)
Staurosporine	< 1.0
Ro 31-8220	< 1.0
H-89	24.8

- **Scientific Information:** P24723 (<http://www.expasy.org/uniprot/P24723>)

