

PKCθ

(Protein kinase C theta type, nPKC-theta, PRKCQ, PRKCT, KPCT)

- **CAT#:** PKCθ
- **Enzyme:** Human PKCθ
- **Substrate:** Histone H1, Lipid activator, 5 μM
ATP 1 μM
- **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	0.8
Ro 31-8220	< 1.0
GW5074	9,900
H-89	5,300

- **Scientific Information:** Q04759 (<http://www.expasy.org/uniprot/Q04759>)

Reference Compound IC50 for PKCθ

