

PKCiota

(Protein kinase C iota type, nPKC-iota, Atypical protein kinase C-lambda/iota, aPKC-lambda/iota, PRKC-lambda/iota, PRKCI, PKCI)

- **CAT#:** PKCiota
- **Enzyme:** Human PKCiota
- **Substrate:** Peptide substrate, [ERM₂PRKRQGSVRRRV], 20 μM
ATP 10 μM
- **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	6.0
Ro 31-8220	94.3
H-89	16,700

- **Scientific Information:** P41743 (<http://www.expasy.org/uniprot/P41743>)

Reference Compound IC50 for PKCiota

