

PKC ζ

(Protein kinase C zeta type, nPKC-zeta, PRKCZ, PKC2, KPCZ)

- **CAT#:** PKC ζ
- **Enzyme:** Human PKC ζ
- **Substrate:** Peptide substrate, [ERM RPRKRQGSVRRRV], 20 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	0.26
Ro 31-8220	1.3
H-89	624

- **Scientific Information:** Q05513 (<http://www.expasy.org/uniprot/Q05513>)

Reference Compound IC50 for PKC ϵ

