

PKC ϵ

(Protein kinase C epsilon type, nPKC-epsilon, PRKCE, PKCE, KPCE)

- **CAT#:** PKC ϵ
- **Enzyme:** Human PKC ϵ
- **Substrate:** Peptide substrate, [ERM ϵ PRKRQGSVRRRV], 20 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 | 0.18 |
| Ro 31-8220 | 0.46 |
| H-89 | 232 |

- **Scientific Information:** Q02156 (<http://www.expasy.org/uniprot/Q02156>)

Reference Compound IC50 for PKC ϵ

