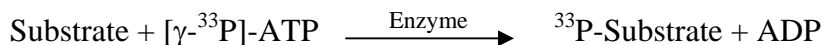


PKCepsilon

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

- **CAT#:** PKCepsilon
- **Enzyme:** Human PKCepsilon
- **Substrate:** Peptide substrate, [ERM_RPRKRQGSVRRRV], 20 μM
ATP 10 μM
- **Reaction:**

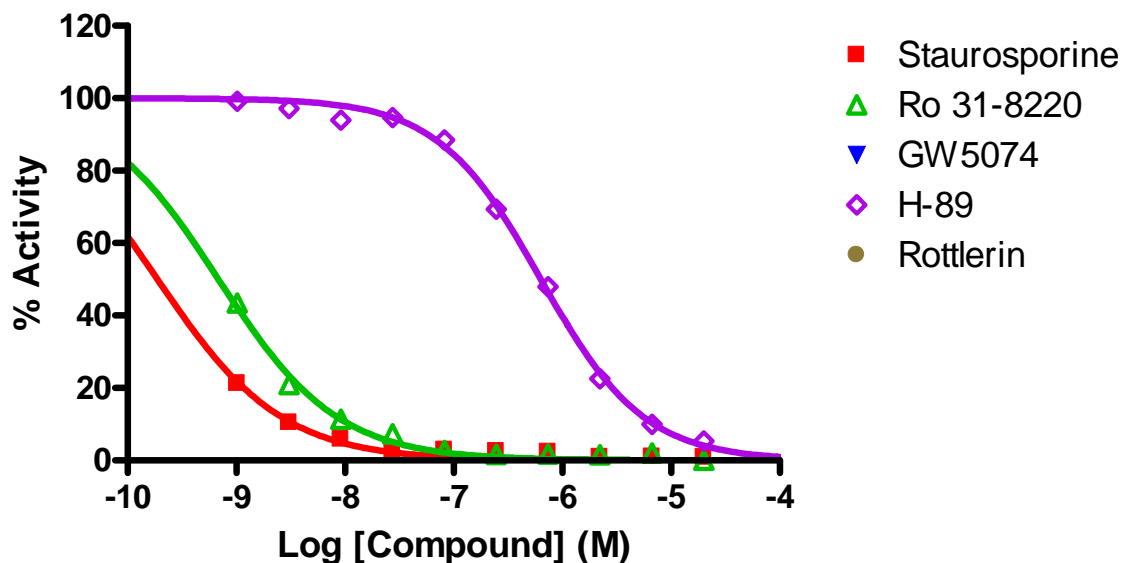


- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 0.19 |
| Ro 31-8220 | 0.69 |
| H-89 | 630 |

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

Reference compound IC50 for PKCepsilon



Service performed
using kinases
under license from  **invitrogen**[™]
part of *life* technologies[™]