

## PKCzeta

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

- **CAT#:** PKCzeta
- **Enzyme:** Human PKCzeta
- **Substrate:** Peptide substrate, [ERM<sub>2</sub>PRKRQGSVRRRV], 20 μM  
ATP 10 μM
- **Reaction:**

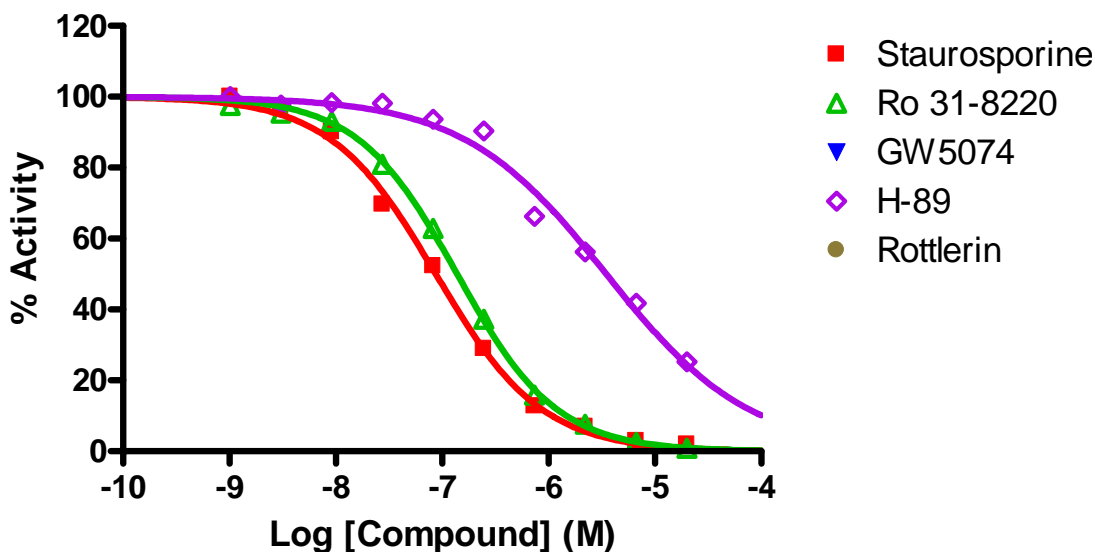


- **Reference Compound IC50s:**

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 87        |
| Ro 31-8220    | 138       |
| H-89          | 3,500     |

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

### Reference compound IC50 for PKCzeta



Service performed  
using kinases  
under license from  **invitrogen**<sup>™</sup>  
part of *life* technologies<sup>™</sup>