

CK1epsilon

(Casein kinase I isoform epsilon, CKI- epsilon, CK1e, CSNK1E, KC1E)

- **CAT#:** CK1epsilon
- **Enzyme:** Human CK1epsilon
- **Substrate:** Peptide substrate, [KRRRAL[pS]VASLPGL], 20 μM
ATP 10 μM
- **Reaction:**

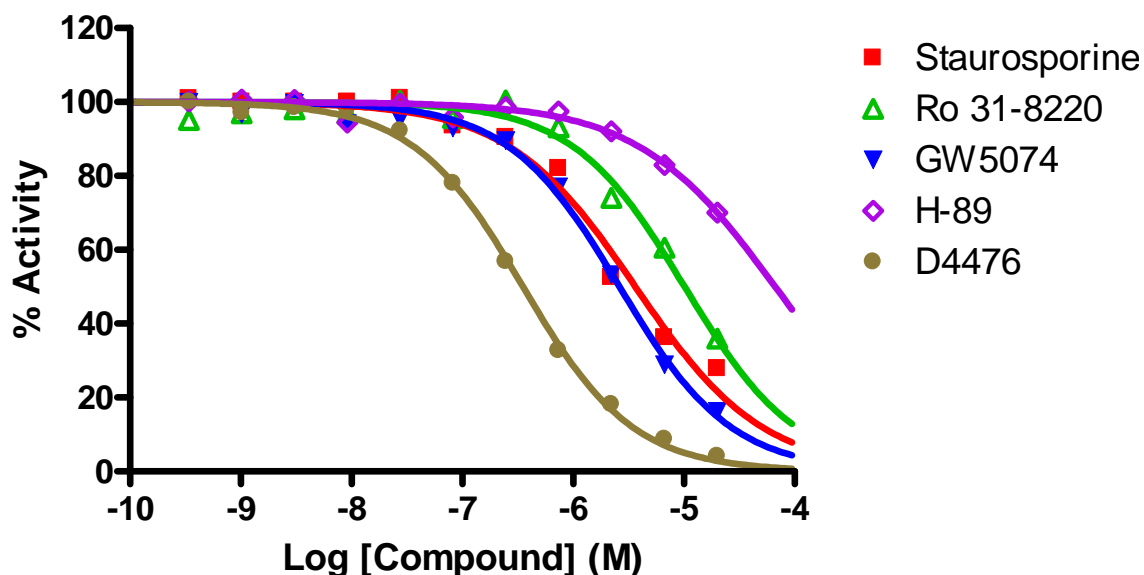


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3,600
Ro 31-8220	10,100
GW 5074	2,600
H-89	66,800
D4476	350

- **Scientific Information:** P49674 (<http://www.expasy.org/uniprot/P49674>)

Reference Compound IC50 for CK1epsilon



Service performed
using kinases
under license from  invitrogen™
part of *Life* technologies™