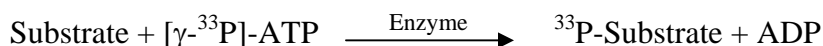


CAMK1a

(Calcium/calmodulin-dependent protein kinase type 1, CaM kinase I, CaM-KI, CaM kinase I alpha, CaMKI-alpha, KCC1A)

- **CAT#:** CAMK1a
- **Enzyme:** Human CAMK1a
- **Substrate:** Peptide substrate, [KKALRRQETVDAL], 20 μ M
Ca²⁺-Calmodulin, 1 μ M
ATP 10 μ M

- **Reaction:**

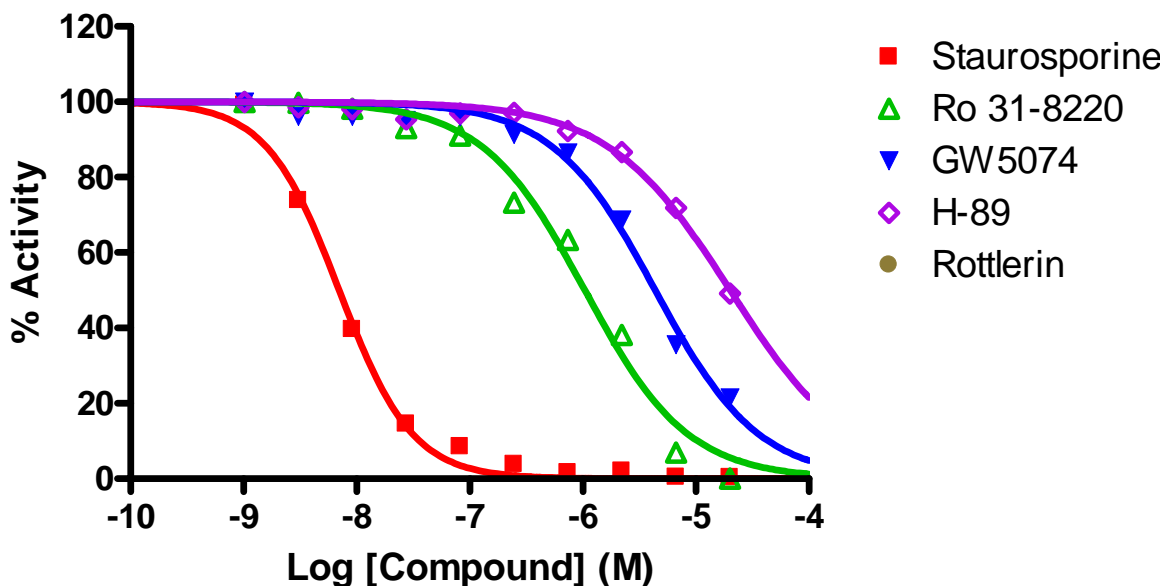


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	7.0
Ro 31-8220	1,000
GW 5074	4,400
H-89	20,000

- **Scientific Information:** Q14012 (<http://www.expasy.org/uniprot/Q14012>)

Reference compound IC50 for CAMK1a



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