

PKCnu/PRKD3

(Protein kinase C nu type, nPKC-nu
Protein kinase EPK2, PRKD3, PRKCN, KPCD3)

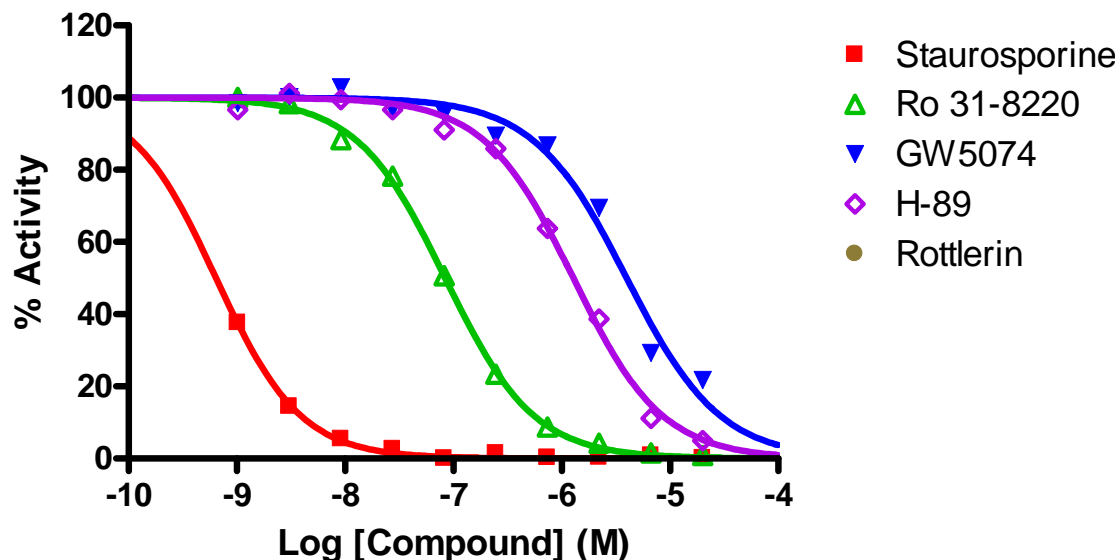
- **CAT#:** PKCnu/PRKD3
- **Enzyme:** Human PKCnu/PRKD3
- **Substrate:** Peptide substrate, [KKLNRTLVA], 20 μ M
ATP 10 μ M
- **Reaction:**

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.6
Ro 31-8220	84
GW 5074	4,000
H-89	1,300

- **Scientific Information:** O94806 (<http://www.expasy.org/uniprot/O94806>)

Reference compound IC50 for PKCnu/PRKD3



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