

## PKG1 $\beta$

(cGMP-dependent protein kinase 1 beta isozyme, cGK 1 beta, cGKI-beta, PRKG1, PRKG1B, PRKGR1B, KGP1B)

- **CAT#:** PKG1 $\beta$
- **Enzyme:** Human PKG1 $\beta$
- **Substrate:** Peptide substrate, [LRRASLG], 100  $\mu$ M  
ATP 1  $\mu$ M
- **Reaction:**  

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

Compounds	IC50 (nM)
Staurosporine	<0.5
Ro 31-8220	25
H-89	6.3
Rottlerin	198,000

- **Scientific Information:** P14619 (<http://www.expasy.org/uniprot/P14619>)

Reference Compound IC50 for PKG1 $\beta$

