

## PAK2

(Serine/threonine-protein kinase PAK 2, p21-activated kinase 2, PAK-2, PAK65, Gamma-PAK, S6/H4 kinase)

- **CAT#:** PAK2
- **Enzyme:** Human PAK2
- **Substrate:** Peptide substrate, [KEAKEKRQEIQIAKRRRLSSLRASTSKSGGSQK], 5  $\mu$ M  
ATP 1  $\mu$ 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	0.35
Ro 31-8220	2,100
GW 5074	19,500
H-89	63,400
Rottlerin	40,000

- **Scientific Information:** Q13177 (<http://www.expasy.org/uniprot/Q13177>)

Reference compound IC50 for PAK2

