

## PAK1

(Serine/threonine-protein kinase PAK 1, p21-activated kinase 1, PAK-1, p65-PAK, Alpha-PAK)

- **CAT#:** PAK1
- **Enzyme:** Human PAK1
- **Substrate:** Peptide substrate, [KEAKEKRQEIQIAKRRRLSSLRASTSKSGGSQK], 20  $\mu$ M  
ATP 10  $\mu$ 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	1.3
Ro 31-8220	7,800
GW 5074	18,600

- **Scientific Information:** Q13153 (<http://www.expasy.org/uniprot/Q13153>)

### Reference compound IC50 for PAK1

