

PAK5

(Serine/threonine-protein kinase PAK 7, p21-activated kinase 7, PAK-7, PAK-5, PAK7, KIAA1264)

- **CAT#:** PAK5
- **Enzyme:** Human PAK5
- **Substrate:** Peptide substrate, [KKLNRTLSFAEPG], 5 μ M
ATP 1 μ 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	2.4
Ro 31-8220	2,700
GW 5074	11,000
H-89	11,300

- **Scientific Information:** Q9P286 (<http://www.expasy.org/uniprot/Q9P286>)

Reference Compound IC50 for PAK5

