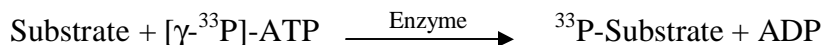


## PAK3

(Serine/threonine-protein kinase PAK 3, p21-activated kinase 3, PAK-3, Beta-PAK, Oligophrenin-3, OPHN3)

- **CAT#:** PAK3
- **Enzyme:** Human PAK3
- **Substrate:** Peptide substrate, [KKLNRTL<sup>S</sup>FAEPG], 5  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.5
Ro 31-8220	898
H-89	10,700

- **Scientific Information:** O75914 (<http://www.expasy.org/uniprot/O75914>)

### Reference Compound IC50 for PAK3

