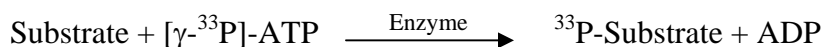


PIM3

(Serine/threonine-protein kinase Pim-3)

- **CAT#:** PIM3
- **Enzyme:** Human PIM3
- **Substrate:** Peptide substrate, [RSRHSSYPAGT], 20 μ M
ATP 1 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.1
Ro 31-8220	9.7
GW 5074	23
H-89	38,300
Rottlerin	2,000

- **Scientific Information:** Q86V86 (<http://www.expasy.org/uniprot/Q86V86>)

Reference compound IC50 for PIM3

