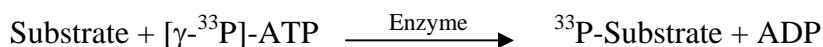


PIM2

(Serine/threonine-protein kinase Pim-2, Pim-2h)

- **CAT#:** PIM2
- **Enzyme:** Human PIM2
- **Substrate:** Peptide substrate, [RSRHSSYPAGT], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	29
Ro 31-8220	750
GW 5074	50
Rottlerin	6,500

- **Scientific Information:** Q9P1W9 (<http://www.expasy.org/uniprot/Q9P1W9>)

Reference compound IC50 for PIM2

