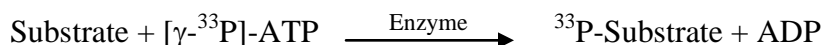


OSR1/OXSR1

(Serine/threonine-protein kinase OSR1, Oxidative stress-responsive 1 protein, KIAA1101)

- **CAT#:** OSR1/OXSR1
- **Enzyme:** Human OSR1/OXSR1
- **Substrate:** Peptide substrate, [RRHYYYDTHNTYYLRTFGHNTRR], 20 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	34
Ro 31-8220	21,000
GW 5074	14,600
Rottlerin	10,200

- **Scientific Information:** O95747 (<http://www.expasy.org/uniprot/O95747>)

Reference compound IC50 for OSR1/OXSR1

