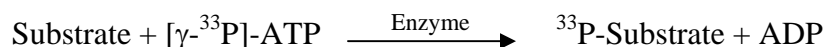


STK38L/NDR2

(Serine/threonine-protein kinase 38-like, NDR2 protein kinase,
Nuclear Dbf2-related kinase 2, KIAA0965, ST38L)

- **CAT#:** STK38L/NDR2
- **Enzyme:** Human STK38L/NDR2
- **Substrate:** Peptide substrate, [KKRNRRLSVA], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.4
Ro 31-8220	1,750
GW 5074	22,000
H-89	810
Rottlerin	>50,000

- **Scientific Information:** Q9Y2H1 (<http://www.expasy.org/uniprot/Q9Y2H1>)

Reference compound IC50 for STK38L/NDR2

