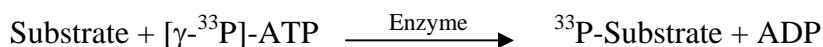


## STK38/NDR1

(Serine/threonine-protein kinase 38, NDR1 protein kinase, Nuclear Dbf2-related kinase 1)

- **CAT#:** STK38/NDR1
- **Enzyme:** Human STK38/NDR1
- **Substrate:** Peptide substrate, [CGRTGRRNSI], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	16
Ro 31-8220	7,200
GW 5074	9,400
H-89	8,000

- **Scientific Information:** Q15208 (<http://www.expasy.org/uniprot/Q15208>)

### Reference compound IC50 for STK38/NDR1

