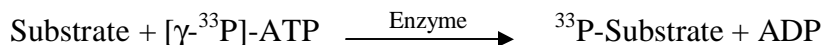


## ROCK2

(Rho-associated protein kinase 2, Rho-associated coiled-coil-containing protein kinase 2, p164 ROCK-2, Rho kinase 2)

- **CAT#:** ROCK2
- **Enzyme:** Human ROCK2
- **Substrate:** Peptide substrate, [KEAKEKRQEIQIAKRRRLSSLRASTSKSGGSQK], 5  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Ro 31-8220	1,400
GW 5074	1,400
H-89	12.5

- **Scientific Information:** O75116 (<http://www.expasy.org/uniprot/O75116>)

### Reference Compound IC50 for ROCK2

