

## STK39/STLK3

(STE20/SPS1-related proline-alanine-rich protein kinase, Ste-20-related kinase, Serine/threonine-protein kinase 39, DCHT, PAK)

- **CAT#:** STK39/STLK3
- **Enzyme:** Human STK39/STLK3
- **Substrate:** Peptide substrate, [RRHYYYDTHNTYLLRTFGHNTRR], 20  $\mu$ M  
ATP 1  $\mu$ M
- **Reaction:**  

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} \text{}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	11.5
Ro 31-8220	21,800
GW 5074	6,900
Rottlerin	8,100

- **Scientific Information:** Q9UEW8 (<http://www.expasy.org/uniprot/Q9UEW8>)

### Reference compound IC50 for STK39/STLK3

