

## TSSK3/STK22C

(Testis-specific serine/threonine-protein kinase 3, TSSK-3, Testis-specific kinase 3, Serine/threonine-protein kinase 22C, SPOGA3, STK22C)

- **CAT#:** TSSK3/STK22C
- **Enzyme:** Human TSSK3/STK22C
- **Substrate:** Peptide substrate, [KKKVSRSGLYRSPSPENLNRPR], 20  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**  
$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	29
Ro 31-8220	14,000
GW 5074	5,500
H-89	>50,000
Rottlerin	>50,000

- **Scientific Information:** Q96PN8 (<http://www.expasy.org/uniprot/Q96PN8>)

### Reference compound IC50 for TSSK3/STK22C

