



# TSSK2,STK22B validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name** : TSSK2,STK22B

**Substrate** : STK S1

**Buffer additives** : 10 mM MgCl<sub>2</sub>, 1 mM DTT

**[TSSK2,STK22B]** : 0.037 ng/μL

**Incubation time at RT** : 50 min

**ATP Km** : 0.2 μM

**Staurosporine IC50** : 157 nM

