



# SIK

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name** : SIK  
**Substrate** : STK S1  
**Buffer additives** : 5 mM MgCl<sub>2</sub>, 1 mM DTT  
**[SIK]** : 1 ng/μL  
**Incubation time at RT** : 60 min  
**ATP Km** : 28 μM  
**Staurosporine IC50** : 3.43 nM

