



# SGK3,SGKL validated with HTRF® KinEASE

## TECHNICAL NOTE

**Enzyme name :** SGK3,SGKL

**Substrate :** STK S3

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

**[SGK3,SGKL] :** 0.001 ng/μL

**Incubation time at RT :** 40 min

**ATP Km :** 13.15 μM

**Staurosporine IC<sub>50</sub> :** 133.5 nM

