



# SGK2

## validated with HTRF® KinEASE

### TECHNICAL NOTE

**Enzyme name :** SGK2

**Substrate :** STK S3

**Buffer additives :** 5 mM MgCl<sub>2</sub>, 1 mM DTT

**[SGK2] :** 0.004 ng/μL

**Incubation time at RT :** 50 min

**ATP Km :** 53.35 μM

**Staurosporine IC<sub>50</sub> :** 56.74 nM

