



SGK1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : SGK1
Substrate : STK S3
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[SGK1] : 0.001 ng/μL
Incubation time at RT : 50 min
ATP Km : 34.54 μM
Staurosporine IC50 : 198.4 nM

