



CHK2,CHEK2 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : CHK2,CHEK2
Substrate : STK S1
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[CHK2,CHEK2] : 0.11 ng/μL
Incubation time at RT : 50 min
ATP Km : n.d. μM
Staurosporine IC50 : 165.1 nM

