



PLK1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PLK1
Substrate : STK S1
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[PLK1] : 1 ng/μL
Incubation time at RT : 60 min
ATP Km : 13.56 μM
Staurosporine IC50 : 17.19 nM

