



Aurora A, AURKA, STK6 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : Aurora A, AURKA, STK6
Substrate : STK S2
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[Aurora A, AURKA, STK6] : 0.064 ng/μL
Incubation time at RT : 40 min
ATP Km : 14.83 μM
Staurosporine IC50 : 19.74 nM

