



CHK1,CHEK1 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : CHK1,CHEK1

Substrate : STK S1

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[CHK1,CHEK1] : 0.01 ng/μL

Incubation time at RT : 50 min

ATP Km : 172 μM

Staurosporine IC₅₀ : 21.43 nM

