



TBK1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : TBK1
Substrate : STK S1
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[TBK1] : 1 ng/μL
Incubation time at RT : 60 min
ATP Km : 18.88 μM
Staurosporine IC₅₀ : 28.62 nM

