



Aurora B, AURKB validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : Aurora B, AURKB

Substrate : STK S2

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[Aurora B, AURKB] : 0.025 ng/μL

Incubation time at RT : 30 min

ATP Km : 1.26 μM

Staurosporine IC₅₀ : 21.73 nM

