



ROCK2,ROK α validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : ROCK2,ROK alpha

Substrate : STK S2

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[ROCK2,ROK α] : 0.16 ng/ μ L

Incubation time at RT : 30 min

ATP Km : 2.75 μ M

Staurosporine IC₅₀ : 4.366 nM

