



PHK γ 2

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PHK gamma2
Substrate : STK S1
Buffer additives : 10 mM MgCl₂, 1 mM DTT
[PHK γ 2] : 0.111 ng/ μ L
Incubation time at RT : 50 min
ATP Km : 2.88 μ M
Staurosporine IC50 : 2.174 nM

