



GPRK4,GRK4 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : GPRK4,GRK4
Substrate : STK S1
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[GPRK4,GRK4] : 1 ng/μL
Incubation time at RT : 90 min
ATP Km : 3.4 μM
Staurosporine IC₅₀ : 25.4 nM

