



GRK6

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : GRK6

Substrate : STK S1

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[GRK6] : 1 ng/μL

Incubation time at RT : 60 min

ATP Km : 9.94 μM

Staurosporine IC₅₀ : 0.68 nM

