



PKAC γ ,PRKACG validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKAC gamma,PRKACG
Substrate : STK S3
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[PKAC γ ,PRKACG] : 0.111 ng/ μ L
Incubation time at RT : 40 min
ATP Km : 5.22 μ M
Staurosporine IC50 : 7.55 nM

