



PKC ι ,PRKCI validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKC ι ,PRKCI

Substrate : STK S1

Buffer additives : 10 mM MgCl₂, 1 mM DTT, lipid activator 1X

[PKC ι ,PRKCI] : 1 ng/ μ L

Incubation time at RT : 20 min

ATP Km : 12.43 μ M

Staurosporine IC₅₀ : 81.79 nM

