



PKC α , PRKCA validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKC alpha, PRKCA

Substrate : STK S1

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

[PKC α , PRKCA] : 0.037 ng/ μ L

Incubation time at RT : 20 min

ATP Km : 2.775 μ M

Staurosporine IC₅₀ : 1.954 nM

