



PKC β 1,PRKCB1 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKC beta1,PRKCB1

Substrate : STK S1

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

[PKC β 1,PRKCB1] : 0.037 ng/ μ L

Incubation time at RT : 25 min

ATP Km : 3.208 μ M

Staurosporine IC50 : 2.525 nM

