



PKC θ , PRKCQ validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKC theta, PRKCQ

Substrate : STK S1

Buffer additives : 10 mM MgCl₂, 1 mM DTT

[PKC θ , PRKCQ] : 0.33 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 7.07 μ M

Staurosporine IC₅₀ : 1.951 nM

