



PKC ϵ ,PRKCE validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKC epsilon,PRKCE

Substrate : STK S1

Buffer additives : 10 mM MgCl₂, 1 mM DTT

[PKC ϵ ,PRKCE] : 0.111 ng/ μ L

Incubation time at RT : 30 min

ATP Km : 5.8 μ M

Staurosporine IC50 : 2.217 nM

