



PKC η , PRKCH validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKC eta, PRKCH

Substrate : STK S1

Buffer additives : 10 mM MgCl₂, 1 mM DTT

[PKC η , PRKCH] : 0.037 ng/ μ L

Incubation time at RT : 50 min

ATP Km : 4.46 μ M

Staurosporine IC50 : 4.663 nM

