



PKC δ , PRKCD validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKC delta, PRKCD

Substrate : STK S1

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

[PKC δ , PRKCD] : 0.111 ng/ μ L

Incubation time at RT : 30 min

ATP Km : 4.059 μ M

Staurosporine IC50 : 2.333 nM

