



PKD2,PRKD2 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PKD2,PRKD2

Substrate : STK S1

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[PKD2,PRKD2] : 0.0014 ng/μL

Incubation time at RT : 50 min

ATP Km : 23.61 μM

Staurosporine IC50 : 9.226 nM

