



PRKX

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PRKX

Substrate : STK S2

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[PRKX] : 0.0123 ng/μL

Incubation time at RT : 30 min

ATP Km : 31.46 μM

Staurosporine IC50 : 6 nM

