



PIM2

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : PIM2
Substrate : STK S3
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[PIM2] : 0.33 ng/μL
Incubation time at RT : 50 min
ATP Km : 3.016 μM
Staurosporine IC50 : 34.5 nM

