



ASK1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : ASK1

Substrate : STK S3

Buffer additives : 5 mM MgCl₂ + 1 mM DTT

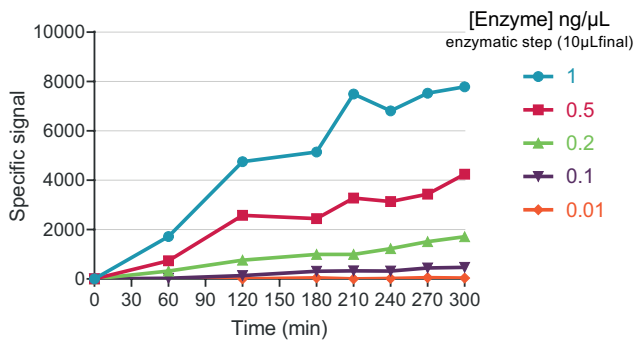
[ASK1] : 1 ng/μL

Incubation time at RT : 210 min

ATP Km : 100 μM

Staurosporine IC50 : 15 nM

ASK1 : Enzyme titration & time course



Staurosporine IC50

