



Haspin,GSG2 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : Haspin,GSG2

Substrate : STK S1

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[Haspin,GSG2] : 1 ng/μL

Incubation time at RT : 90 min

ATP Km : 7.6 μM

Staurosporine IC50 : 0.84 nM

