



YES1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : YES1

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[YES1] : 0.111 ng/μL

Incubation time at RT : 30 min

ATP Km : 24.38 μM

Staurosporine IC50 : 27.58 nM

