



TIE2,TEK validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : TIE2,TEK

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM MnCl₂, 1 mM DTT, SEB

[TIE2,TEK] : 0.111 ng/μL

Incubation time at RT : 50 min

ATP Km : 6.4 μM

Staurosporine IC50 : 93 nM

