



SRC

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : SRC
Substrate : TK
Buffer additives : 5 mM MgCl₂, 1 mM DTT
[SRC] : 0.111 ng/μL
Incubation time at RT : 50 min
ATP Km : 19.02 μM
Staurosporine IC₅₀ : 52.9 nM

