



SYK

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : SYK

Substrate : TK

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

[SYK] : 0.0123 ng/μL

Incubation time at RT : 30 min

ATP Km : 0.11 μM

Staurosporine IC₅₀ : 2.66 nM

