



RET

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : RET

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[RET] : 0.037 ng/μL

Incubation time at RT : 30 min

ATP Km : 7.915 μM

Staurosporine IC₅₀ : 675 nM

