



FGFR1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FGFR1

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[FGFR1] : 0.0123 ng/μL

Incubation time at RT : 50 min

ATP Km : 99.99 μM

Staurosporine IC₅₀ : 37.89 nM

