



FGFR3

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FGFR3

Substrate : TK

Buffer additives : 5 mM MgCl₂, 1 mM DTT

[FGFR3] : 0.0123 ng/μL

Incubation time at RT : 30 min

ATP Km : 38.88 μM

Staurosporine IC₅₀ : 32.2 nM

