



FLT4, VEGFR3 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FLT4, VEGFR3

Substrate : TK

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

[FLT4, VEGFR3] : 0.0123 ng/μL

Incubation time at RT : 30 min

ATP Km : 0.4 μM

Staurosporine IC₅₀ : 4.389 nM

