



FLT1, VEGFR1 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FLT1, VEGFR1

Substrate : TK

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

[FLT1, VEGFR1] : 0.037 ng/μL

Incubation time at RT : 50 min

ATP Km : 3.88 μM

Staurosporine IC₅₀ : 11.97 nM

