



EPHB1

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : EPHB1

Substrate : TK

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

[EPHB1] : 0.037 ng/μL

Incubation time at RT : 40 min

ATP Km : 0.69 μM

Staurosporine IC₅₀ : 100.4 nM

