



EPHA7

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : EPHA7

Substrate : TK

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

[EPHA7] : 0.037 ng/μL

Incubation time at RT : 30 min

ATP Km : 0.503 μM

Staurosporine IC50 : 99.27 nM

