



EPHA5 validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : EPHA5

Substrate : TK

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

[EPHA5] : 0.037 ng/μL

Incubation time at RT : 50 min

ATP Km : 4.8 μM

Staurosporine IC50 : 3230 nM

