



EPHA3

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : EPHA3

Substrate : TK

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

[EPHA3] : 0.111 ng/μL

Incubation time at RT : 50 min

ATP Km : 7.9 μM

Staurosporine IC₅₀ : 3242 nM

