



FER

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FER

Substrate : TK

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

[FER] : 0.333 ng/μL

Incubation time at RT : 50 min

ATP Km : 5.3 μM

Staurosporine IC50 : 168 nM

