



FGR

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : FGR

Substrate : TK

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

[FGR] : 0.01 ng/μL

Incubation time at RT : 50 min

ATP Km : 8.8 μM

Staurosporine IC50 : 135.2 nM

