



ABL2,Arg validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : ABL2,Arg

Substrate : TK

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

[ABL2,Arg] : 0.111 ng/μL

Incubation time at RT : 50 min

ATP Km : 25.9 μM

Staurosporine IC50 : 1727 nM

