



BTK

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : BTK

Substrate : TK

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

[BTK] : 0.111 ng/μL

Incubation time at RT : 50 min

ATP Km : 28.2 μM

Staurosporine IC50 : 2640 nM

