



BMX

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : BMX

Substrate : TK

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

[BMX] : 0.333 ng/μL

Incubation time at RT : 50 min

ATP Km : 26.4 μM

Staurosporine IC50 : 1731 nM

