



AXL

validated with HTRF® KinEASE

TECHNICAL NOTE

Enzyme name : AXL

Substrate : TK

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

[AXL] : 0.111 ng/μL

Incubation time at RT : 50 min

ATP Km : 21 μM

Staurosporine IC50 : 252.5 nM

