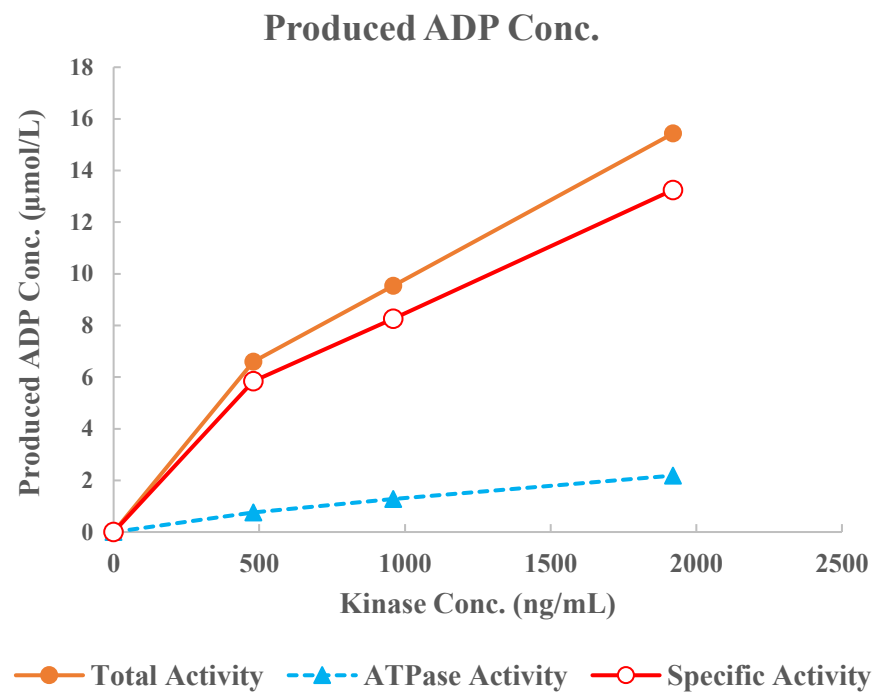
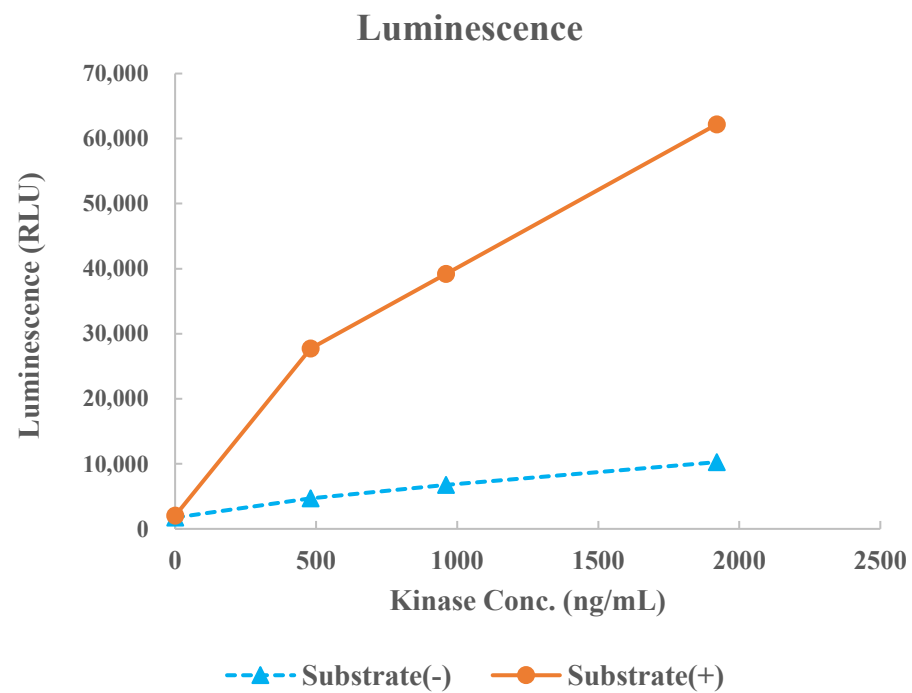


MST1(STK4)



Kinase	MST1(STK4), 07-116
Substrate	Biotin-IRS1, BTSU07
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	20 μmol/L Cantharidin
Reaction Time	1 hour