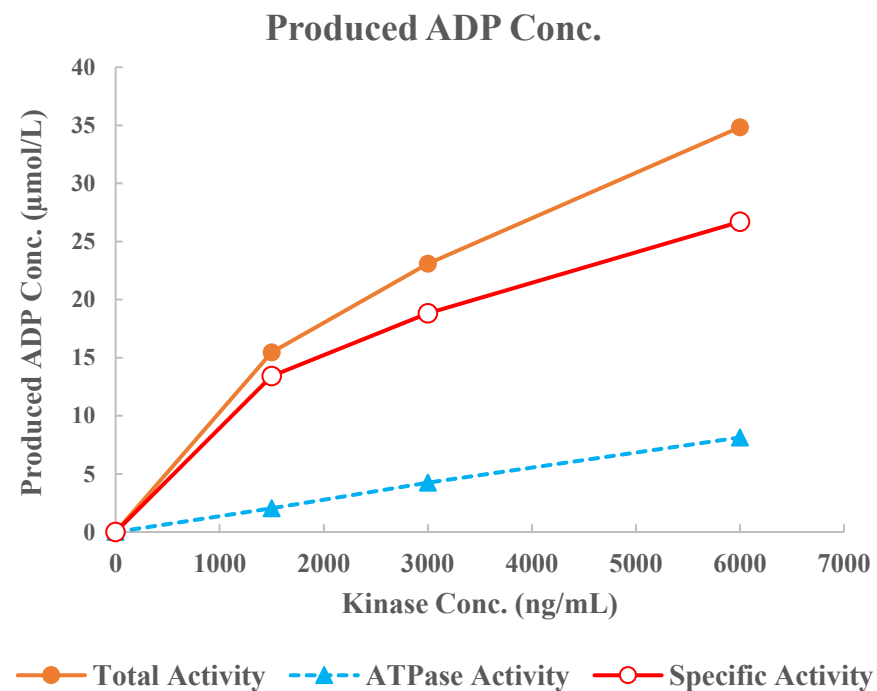
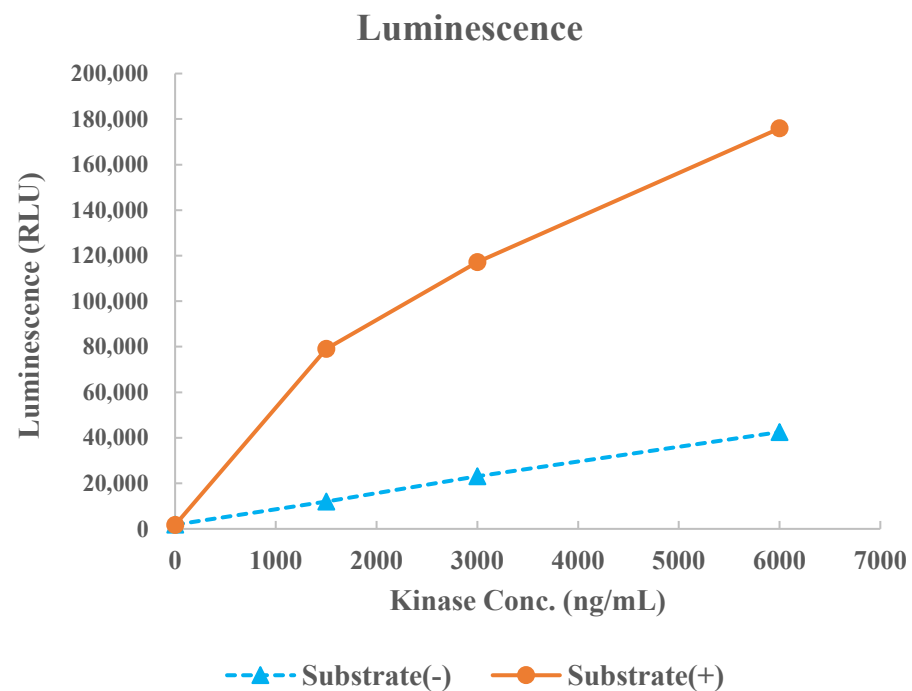


MST2(STK3)



Kinase	MST2(STK3), 07-117
Substrate	Biotin-IRS1, BTSU07
Substrate Conc.	50 μmol/L
ATP Conc.	75 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	10 μmol/L Cantharidin
Reaction Time	1 hour