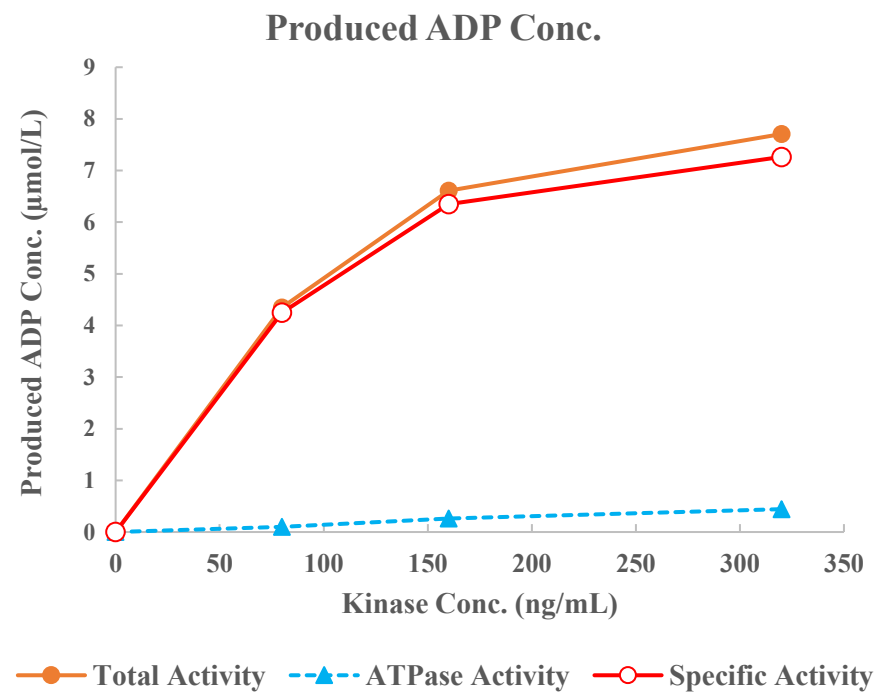
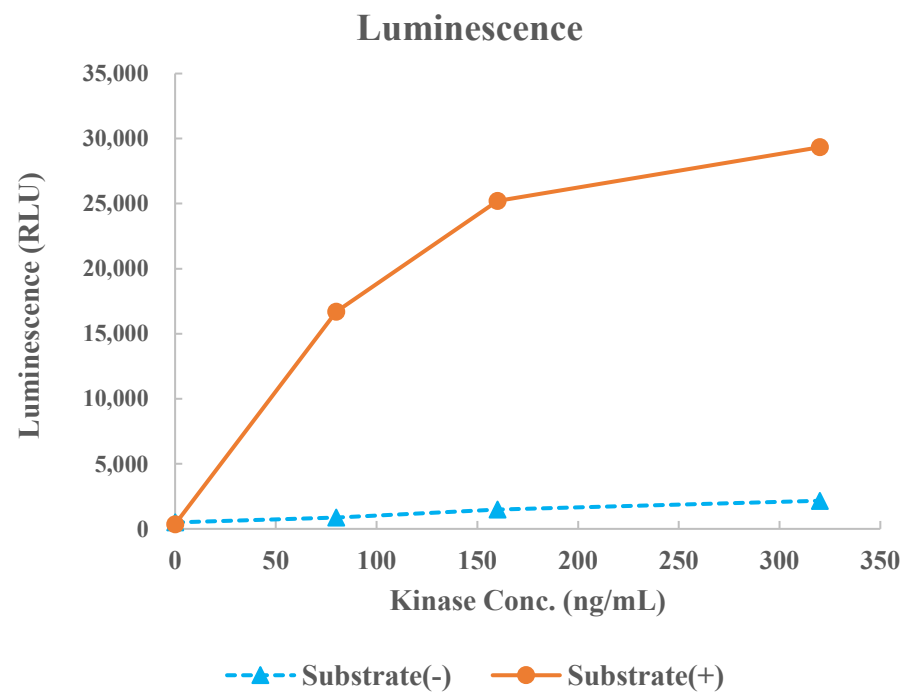


SRPK2



Kinase	SRPK2, 04-161
Substrate	Biotin-DYRKtide, BTSU06
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour