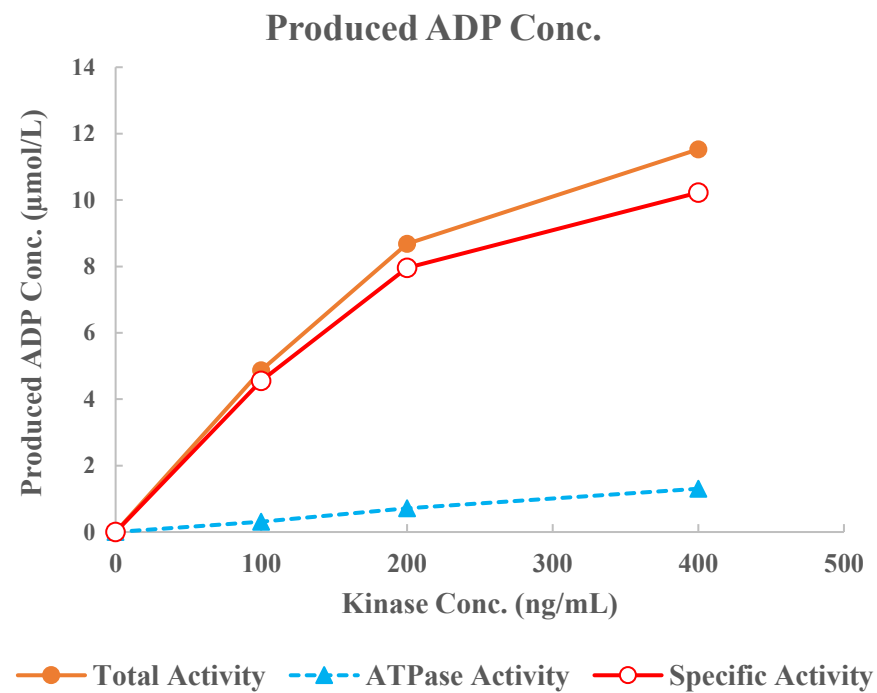
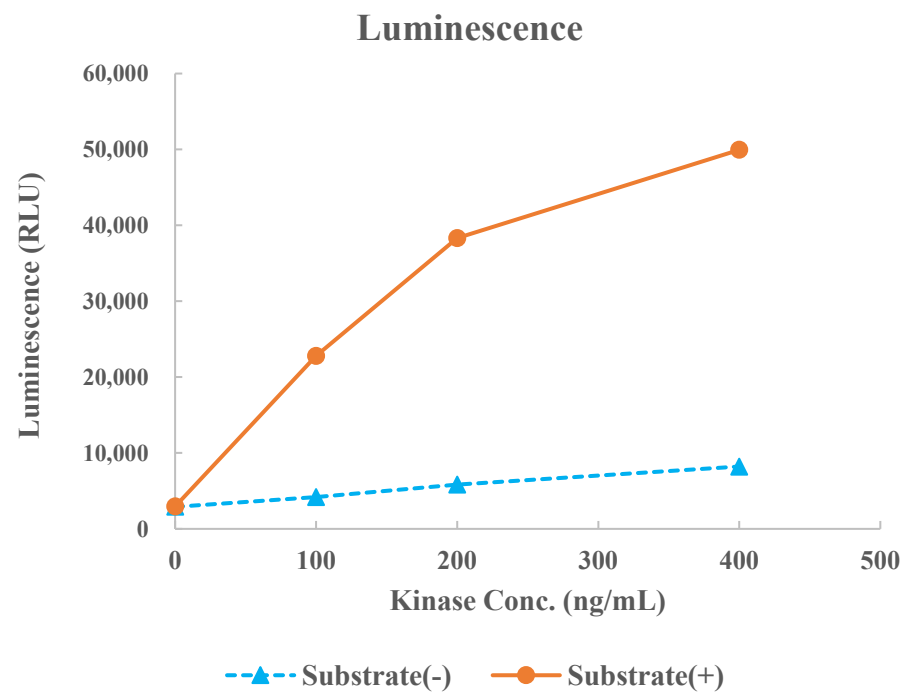


CLK2



Kinase	CLK2, 04-127
Substrate	Biotin-DYRKtide, BTSU06
Substrate Conc.	10 μmol/L
ATP Conc.	150 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour