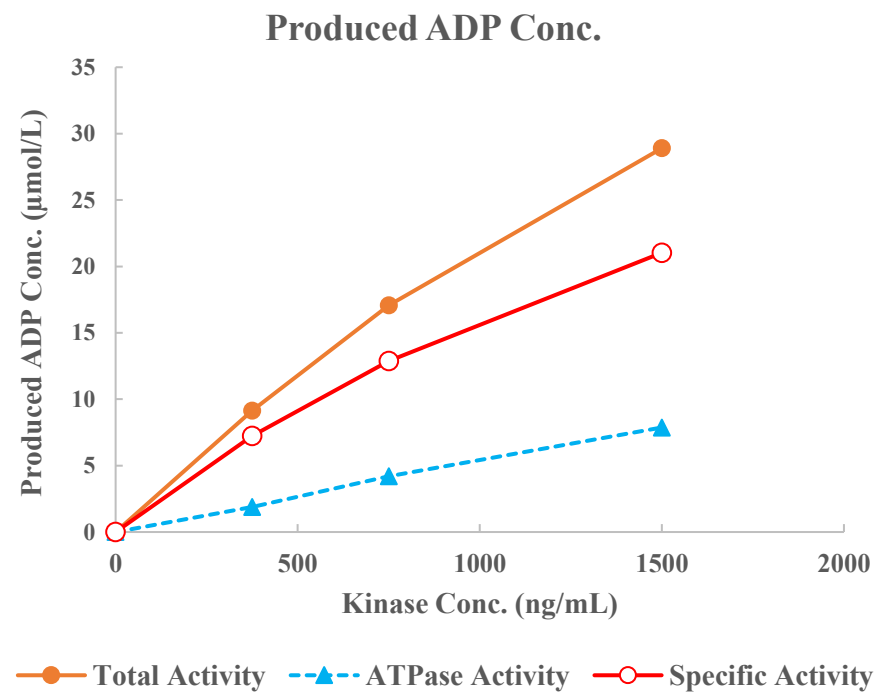
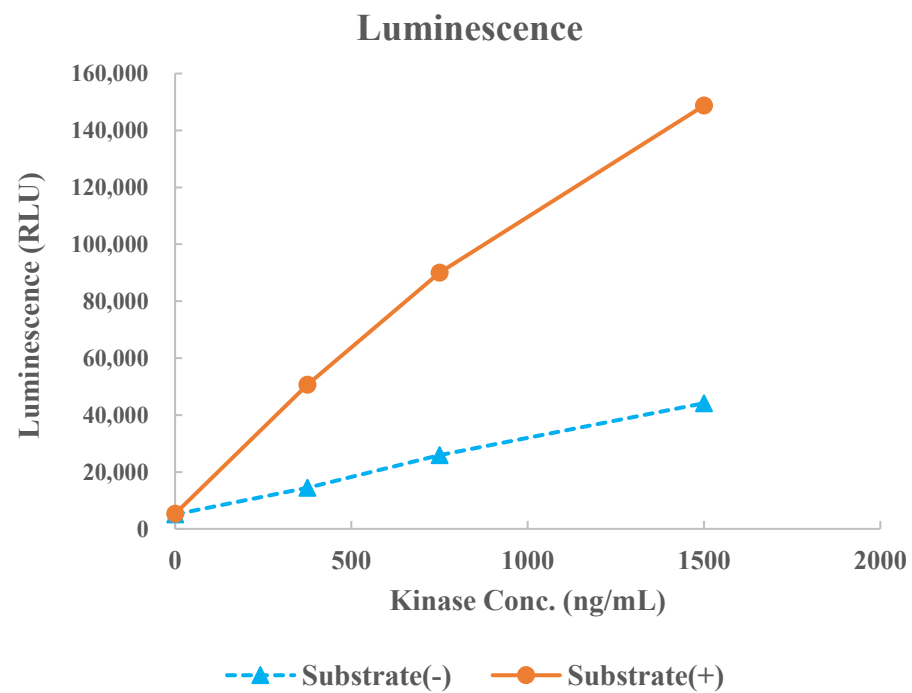


CLK3



Kinase	CLK3, 04-128
Substrate	Biotin-DYRKtide, BTSU06
Substrate Conc.	50 μmol/L
ATP Conc.	75 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour