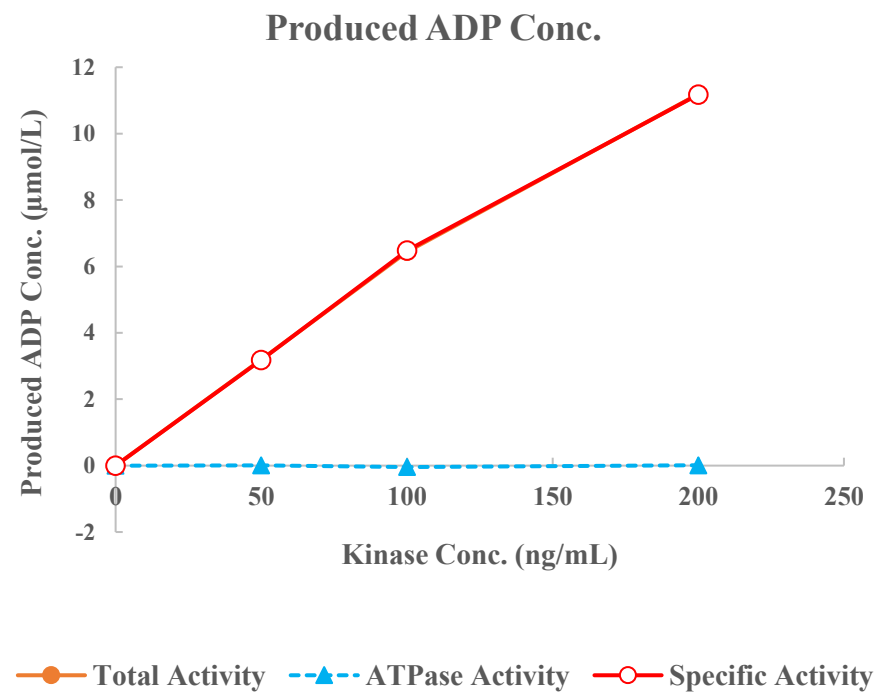
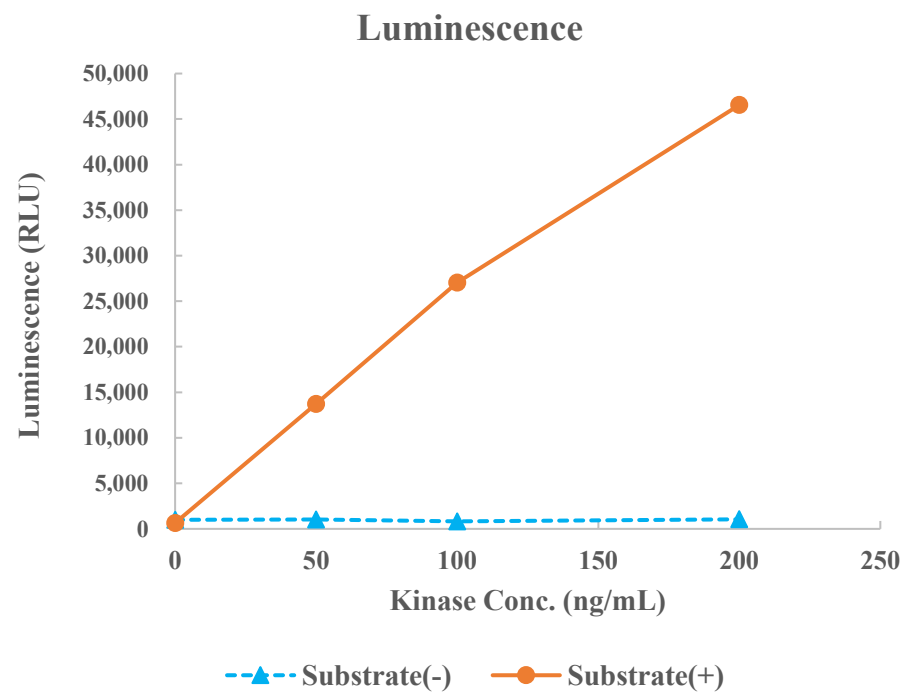


CGK2(PRKG2)



Kinase	CGK2(PRKG2), 01-143
Substrate	Biotin-Kemptide, BTSU09
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	5 μmol/L cGMP
Reaction Time	1 hour