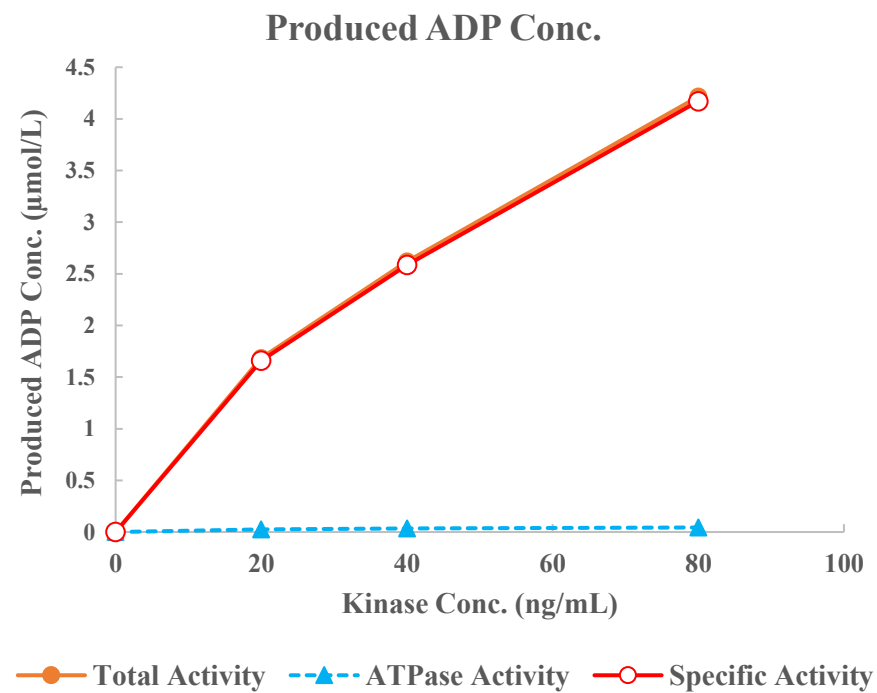
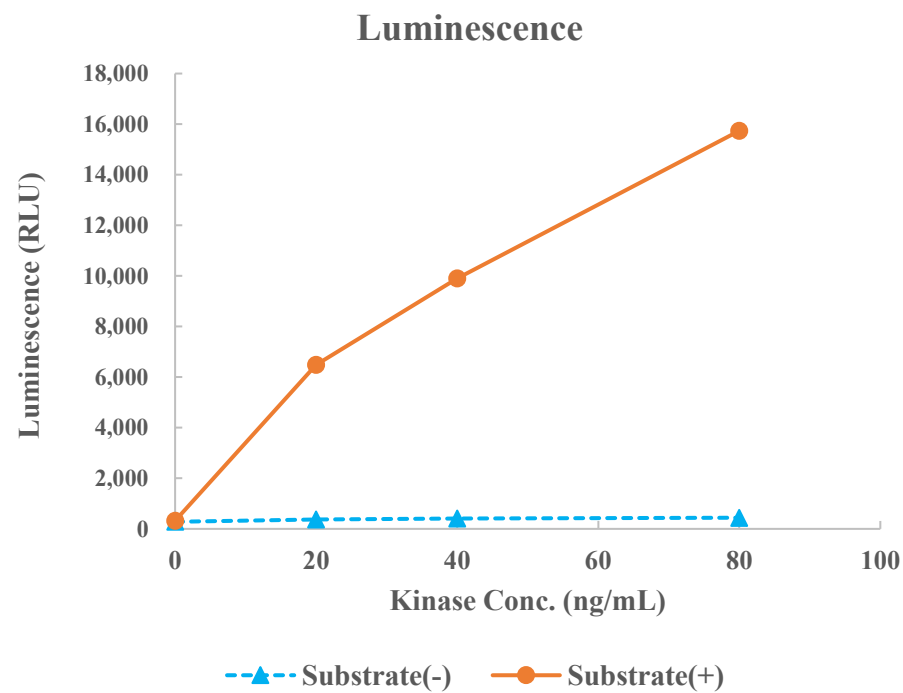


PKAC β (PRKACB)



Kinase	PKAC β (PRKACB), 01-128
Substrate	Biotin-Kemptide, BTSU09
Substrate Conc.	10 μ mol/L
ATP Conc.	5 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour