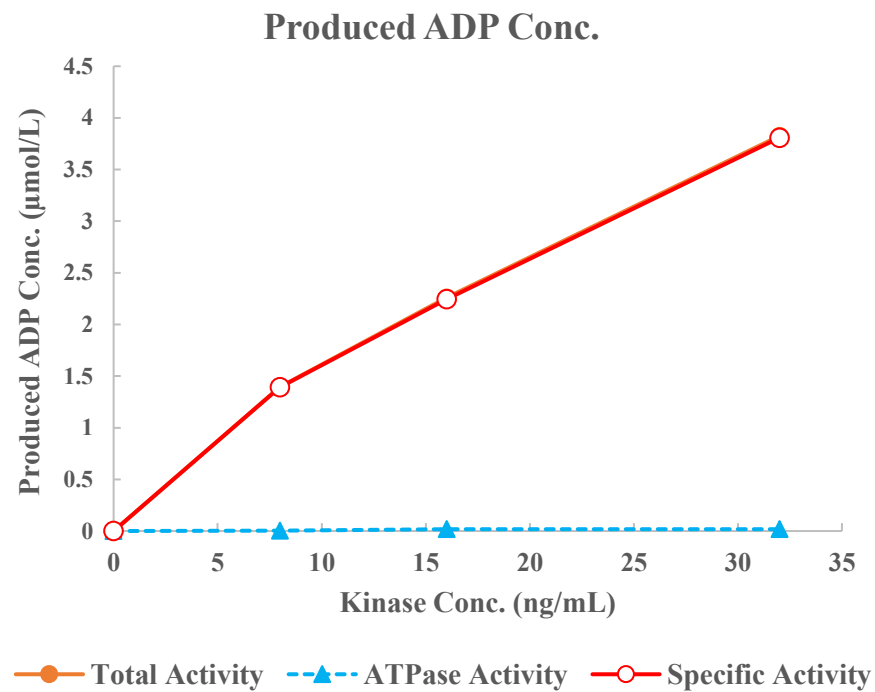
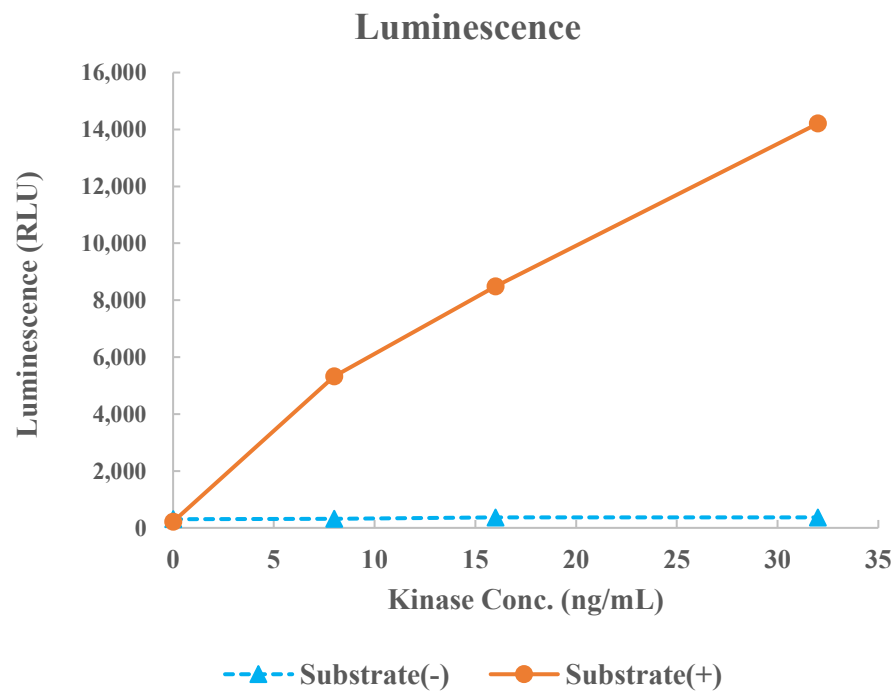


PKAC α (PRKACA)



Kinase	PKAC α (PRKACA), 01-127
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