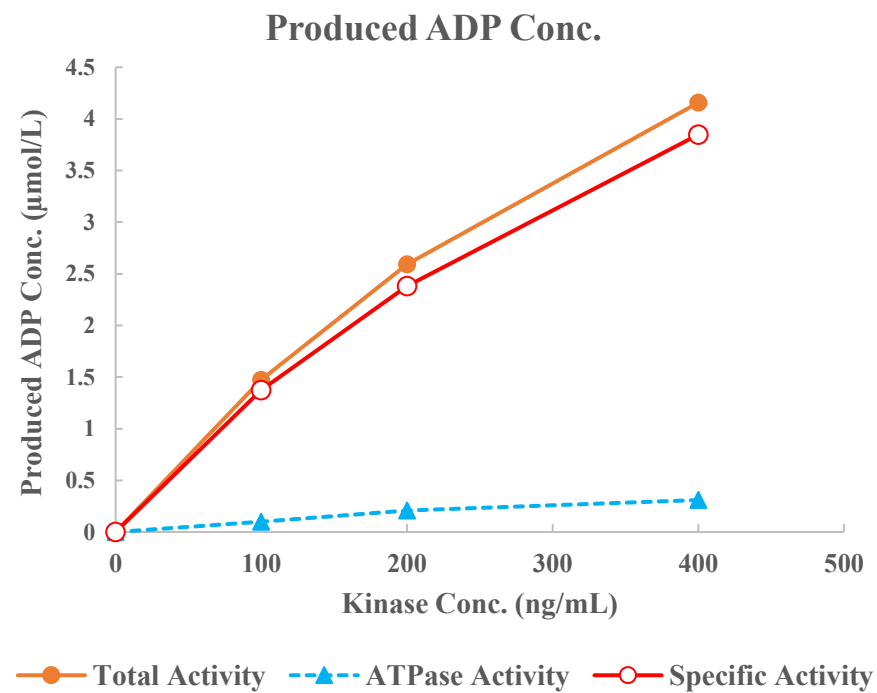
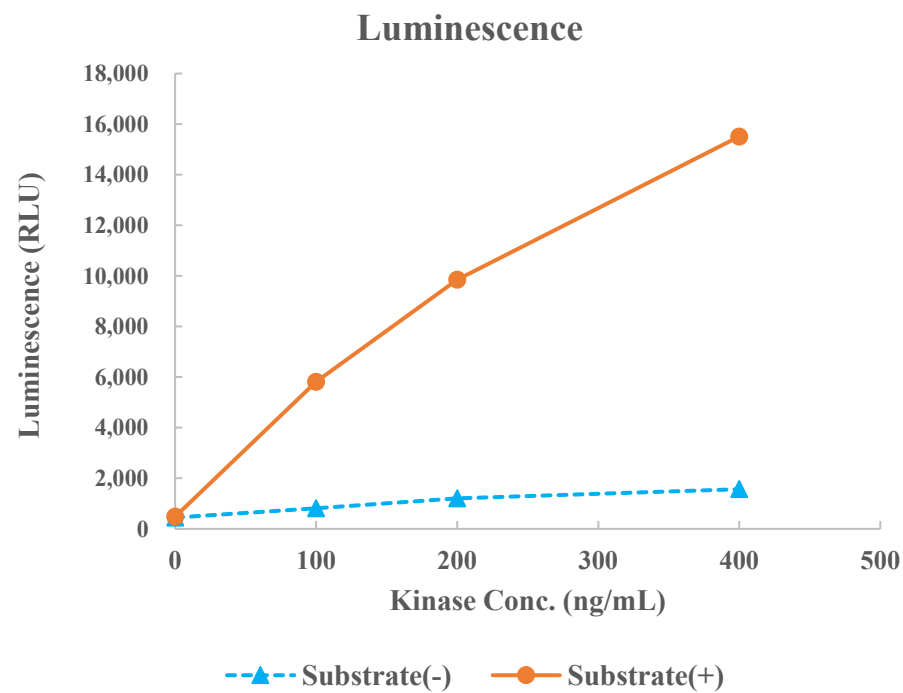


GSK3 α (GSK3A)



Kinase	GSK3 α (GSK3A), 04-140
Substrate	Biotin-CREBtide-p, BTSU22
Substrate Conc.	10 μ mol/L
ATP Conc.	10 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour