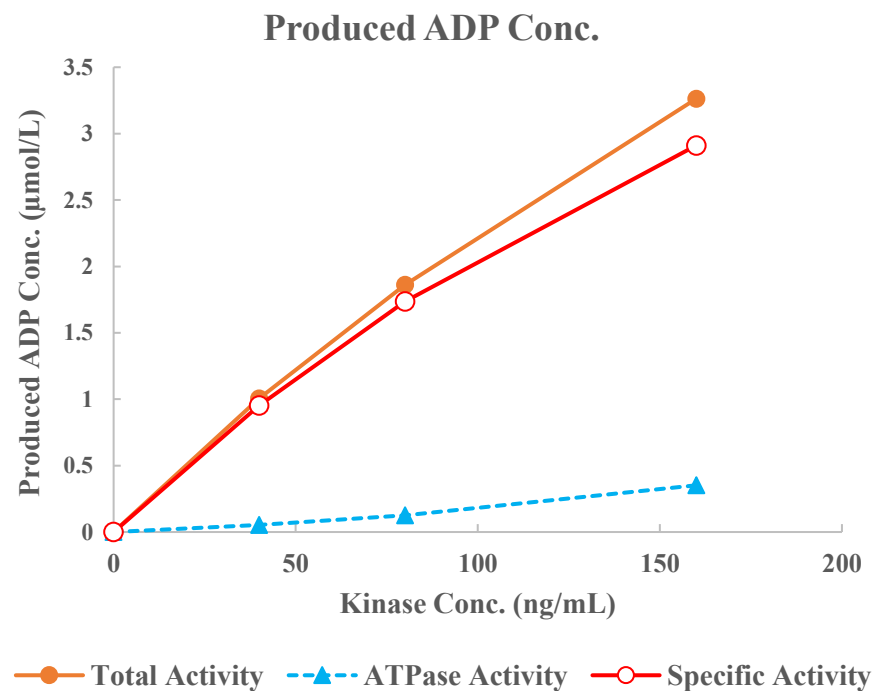
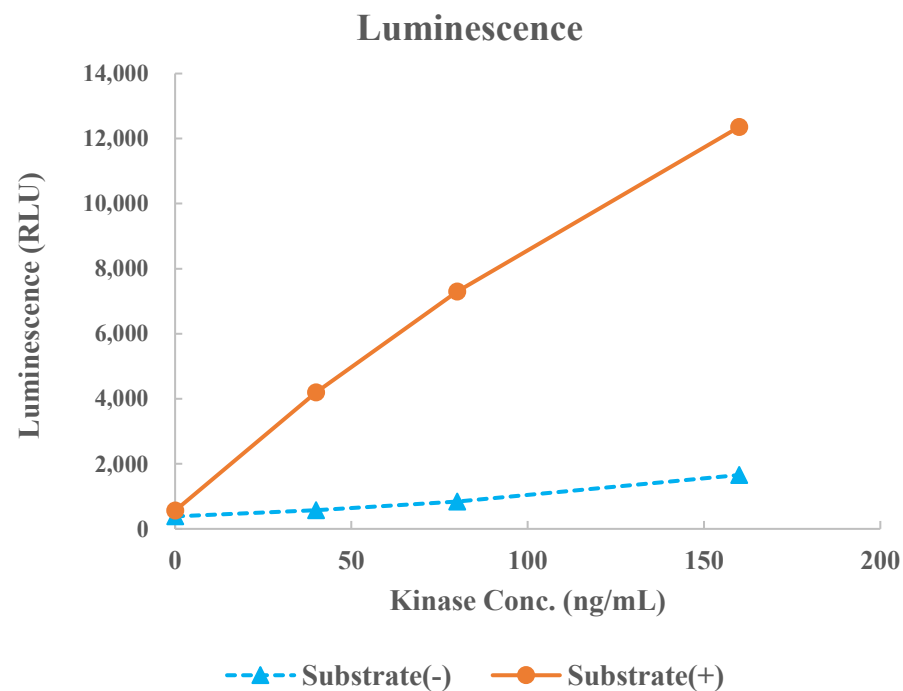


GSK3 β (GSK3B)



Kinase	GSK3 β (GSK3B), 04-141
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