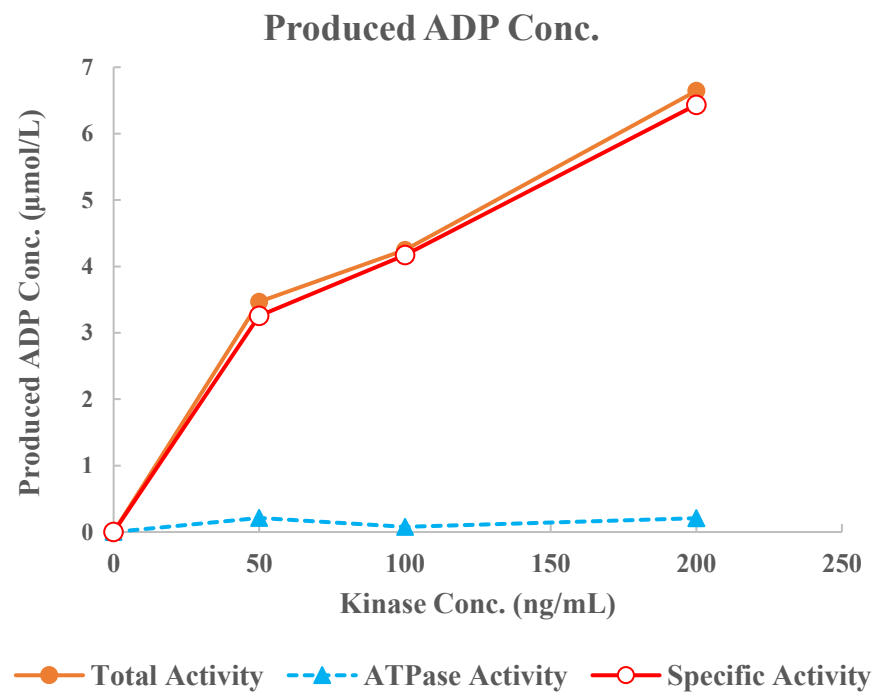
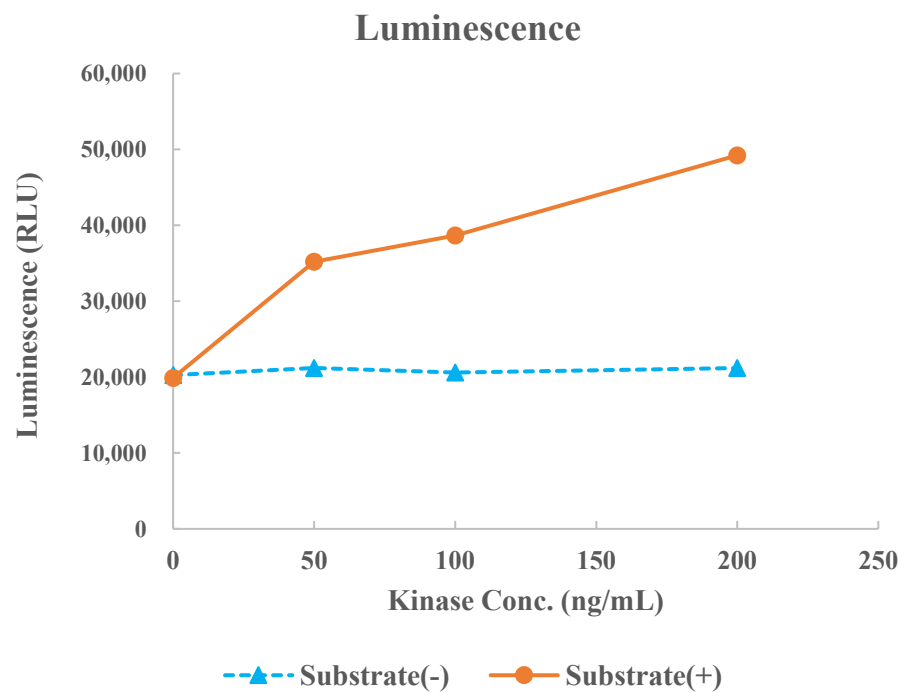


CaMK1 α (CAMK1)



Kinase	CaMK1 α (CAMK1), 02-104
Substrate	Biotin-GS peptide, BTSU05
Substrate Conc.	10 μ mol/L
ATP Conc.	1000 μ mol/L
Metal	Mg, Ca
Metal Conc.	5, 1 mmol/L
Additives	10 μ g/mL Calmodulin
Reaction Time	1 hour