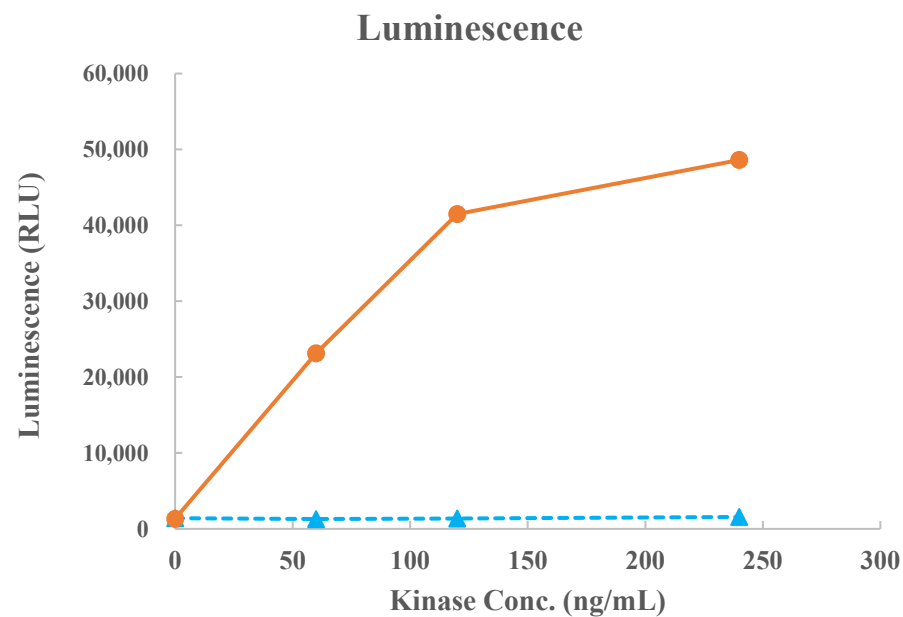
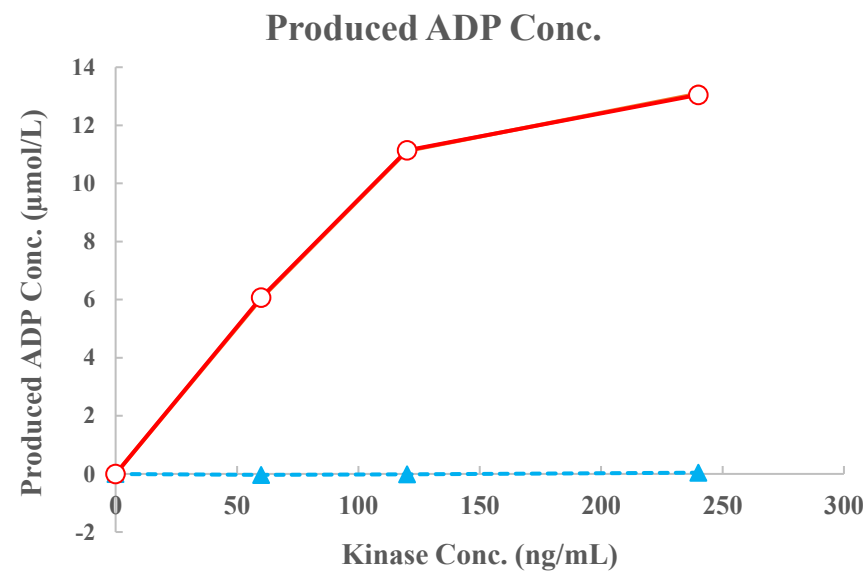


CaMK2 α (CAMK2A)



--▲-- Substrate(-) ● Substrate(+)



● Total Activity --▲-- ATPase Activity ○ Specific Activity

Kinase	CaMK2 α (CAMK2A), 02-109
Substrate	Biotin-GS peptide, BTSU05
Substrate Conc.	10 μ mol/L
ATP Conc.	50 μ mol/L
Metal	Mg, Ca
Metal Conc.	5, 1 mmol/L
Additives	10 μ g/mL Calmodulin
Reaction Time	1 hour