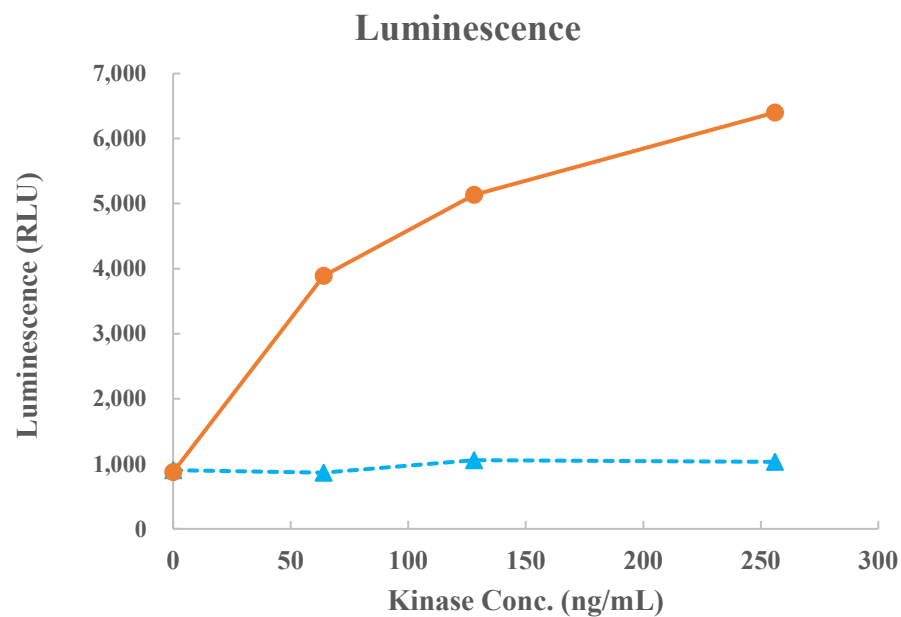
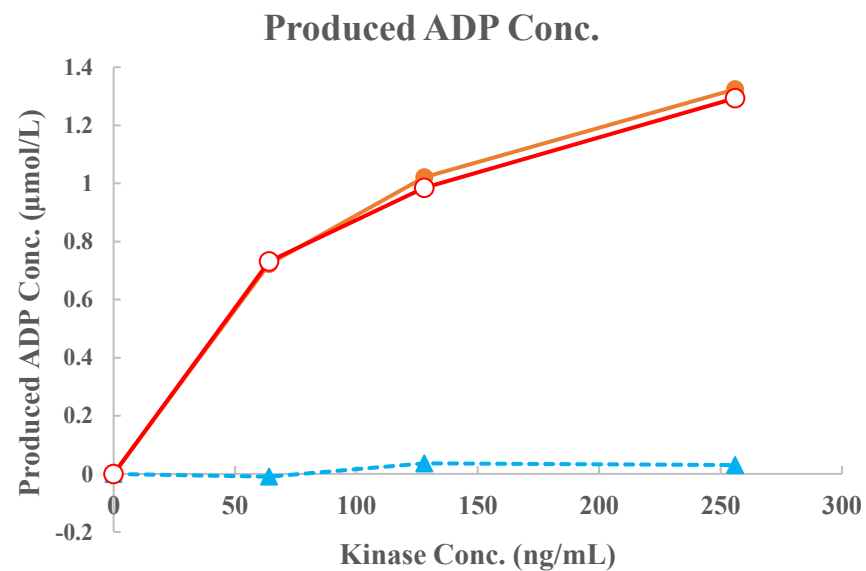


PKC α (PRKCI)



--▲-- Substrate(-) ● Substrate(+)



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

Kinase	PKC α (PRKCI), 01-139
Substrate	Biotin-PKC peptide, BTSU10
Substrate Conc.	10 μ mol/L
ATP Conc.	25 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour