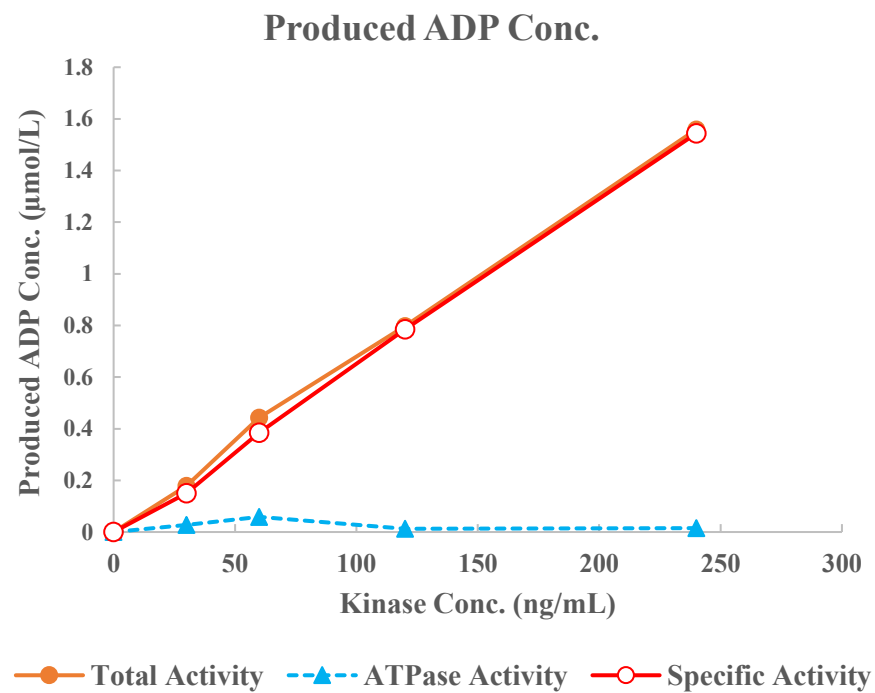
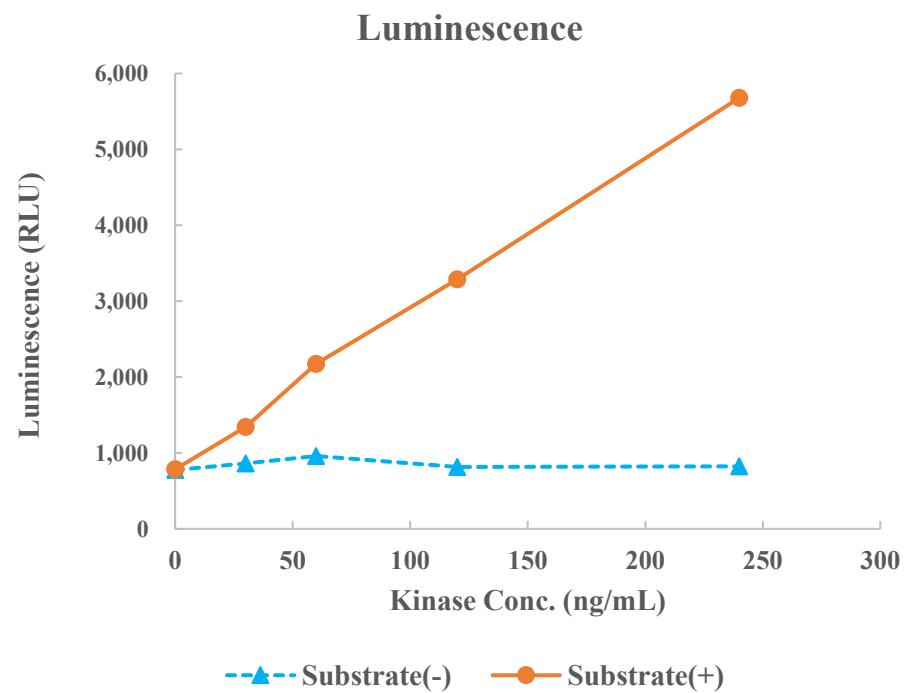


PKC ζ (PRKCZ)



Kinase	PKC ζ (PRKCZ), 01-141
Substrate	Biotin-PKC peptide, BTSU10
Substrate Conc.	10 μ mol/L
ATP Conc.	10 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour