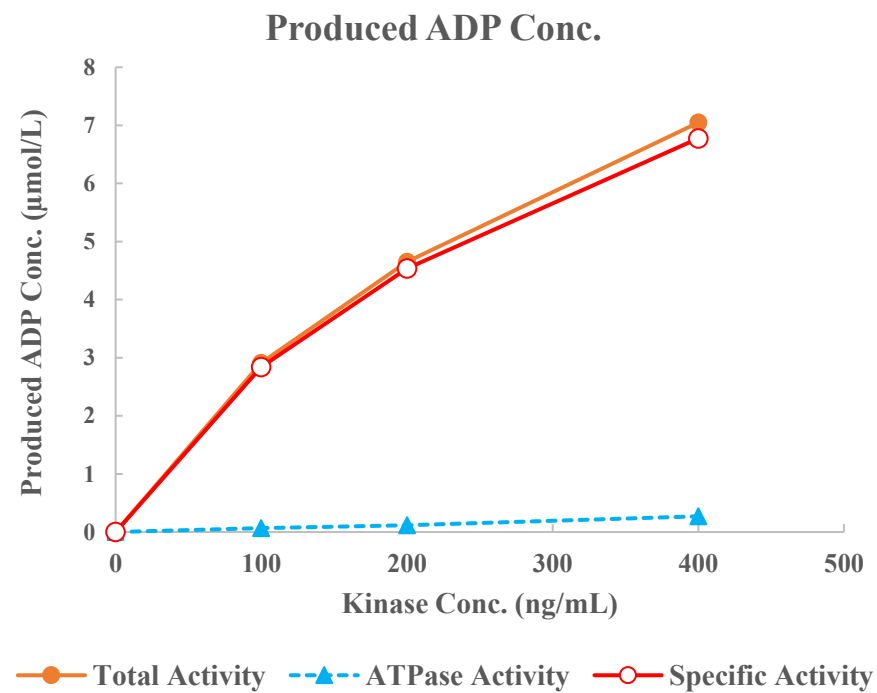
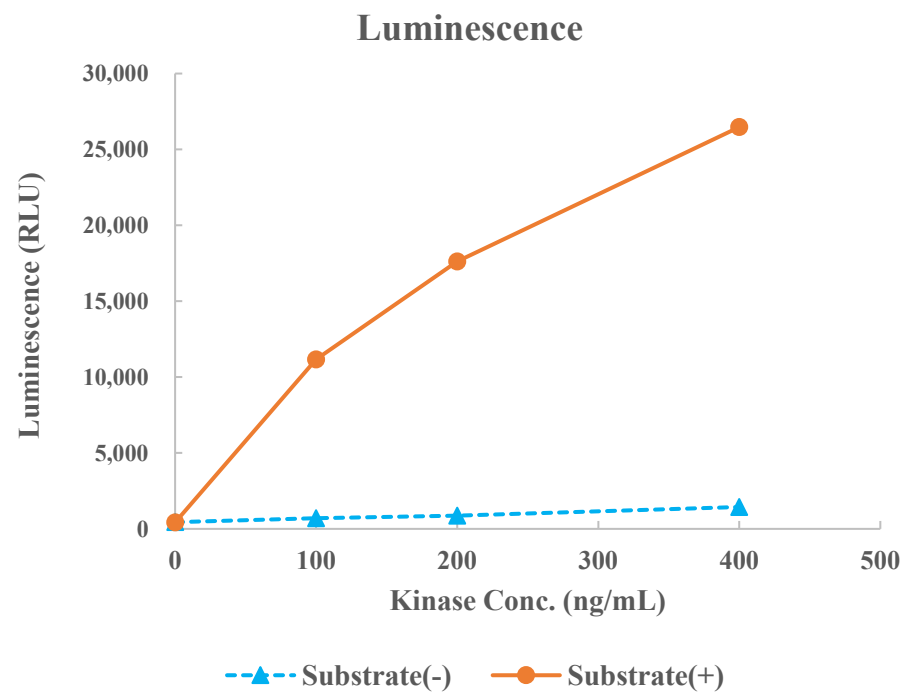


CK1 γ 2(CSNK1G2)



Kinase	CK1 γ 2(CSNK1G2), 03-106
Substrate	Biotin-CKtide, BTSU12
Substrate Conc.	10 μ mol/L
ATP Conc.	10 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour