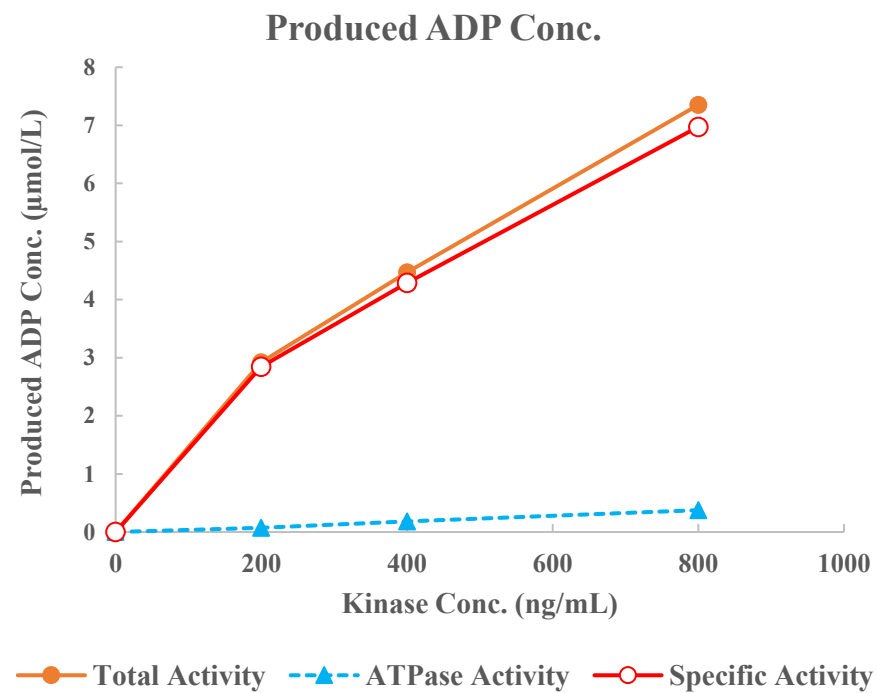
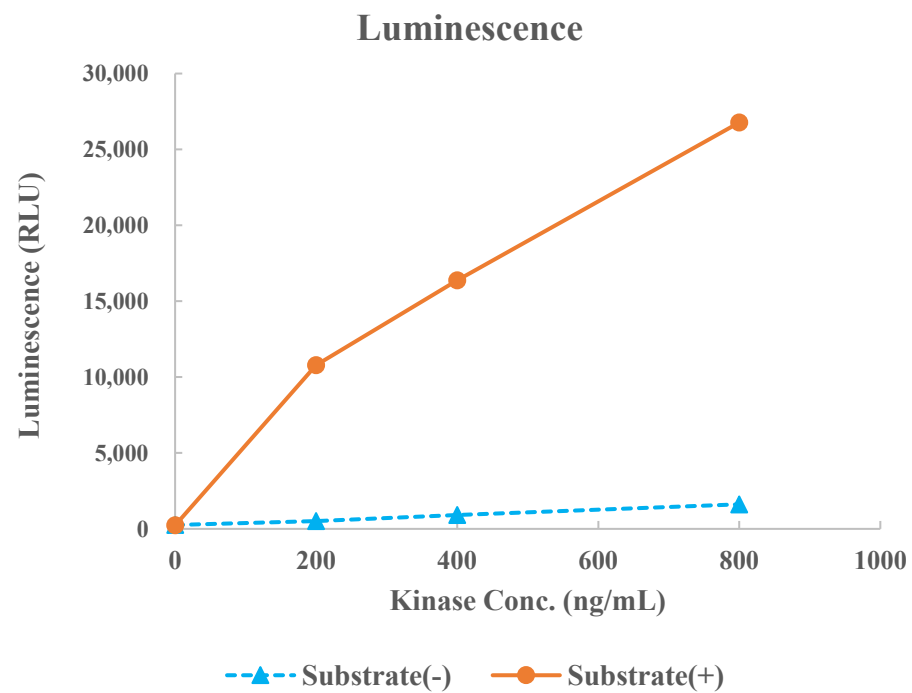


CK1 γ 3(CSNK1G3)



Kinase	CK1 γ 3(CSNK1G3), 03-107
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