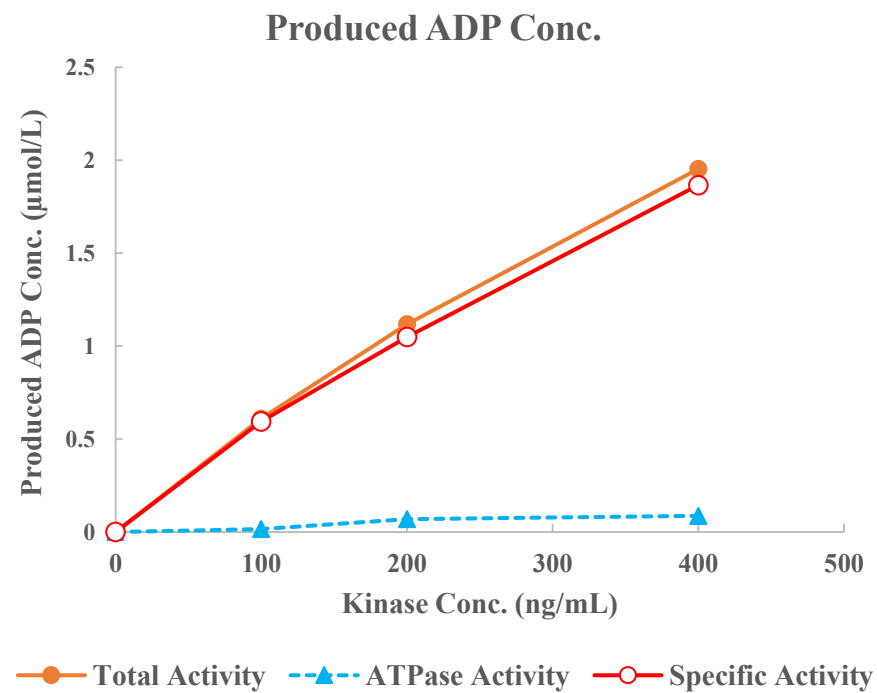
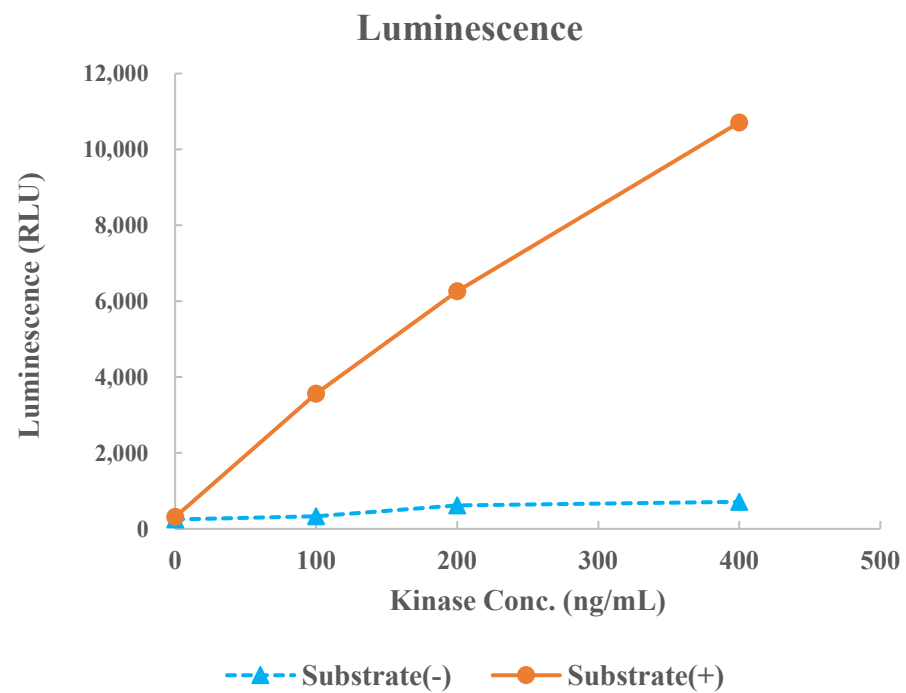


CK1 α (CSNK1A1)



Kinase	CK1 α (CSNK1A1), 03-101
Substrate	Biotin-CKtide, BTSU12
Substrate Conc.	50 μ mol/L
ATP Conc.	5 μ mol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour