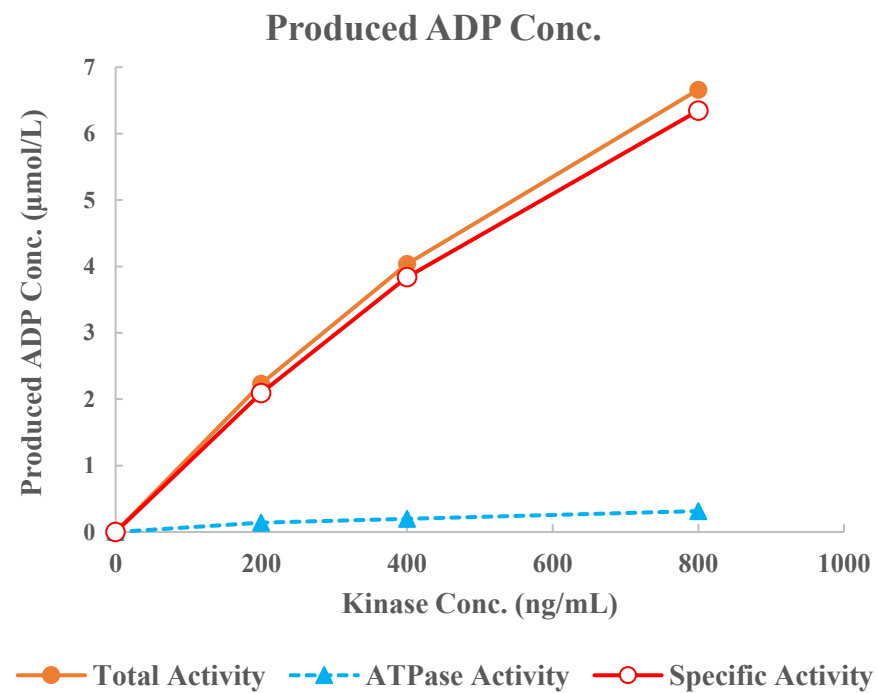
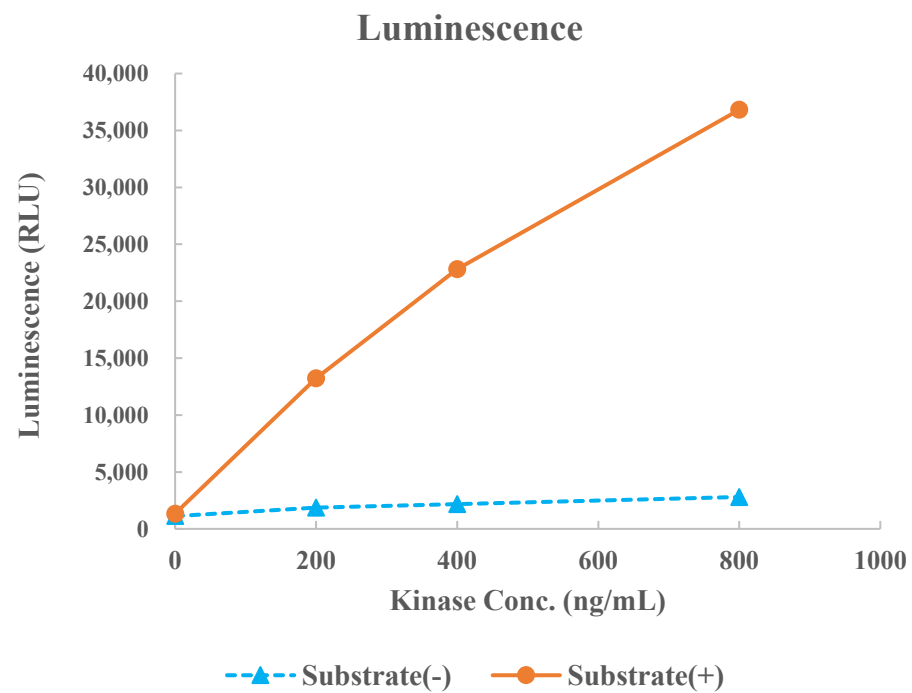


CK1ε(CSNK1E)



Kinase	CK1ε(CSNK1E), 03-104
Substrate	Biotin-CKtide, BTSU12
Substrate Conc.	50 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour