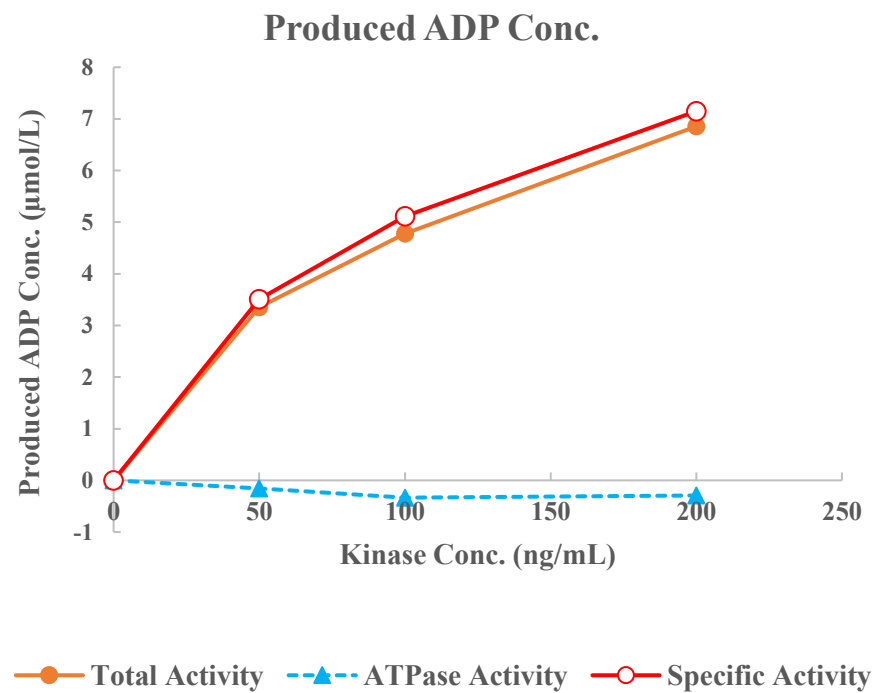
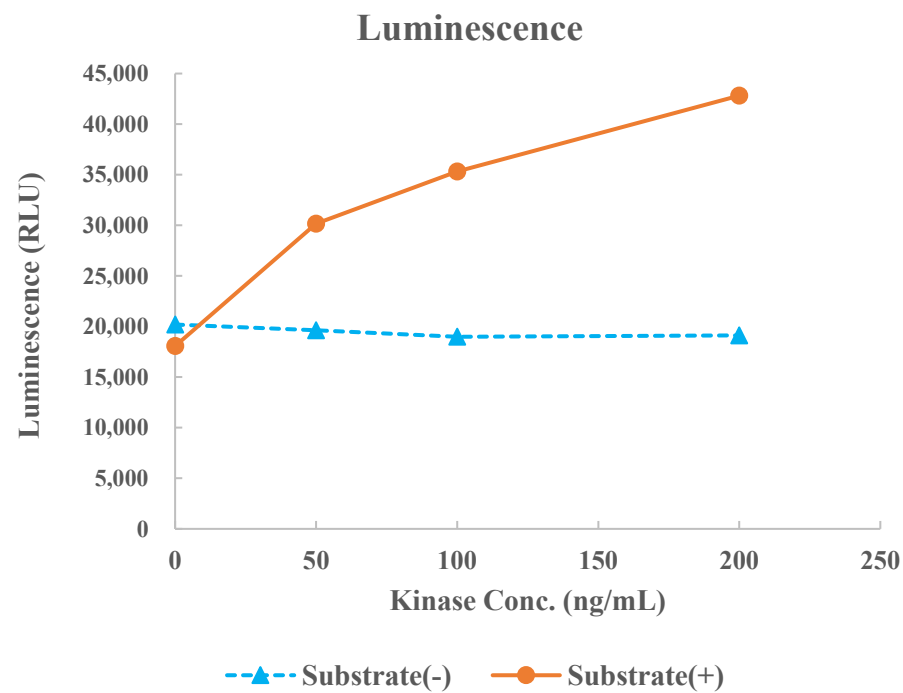


CDK3/CycE1



Kinase	CDK3/CycE1, 04-104
Substrate	Biotin-Modified Histone H1, BTSU13
Substrate Conc.	10 μmol/L
ATP Conc.	1000 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour