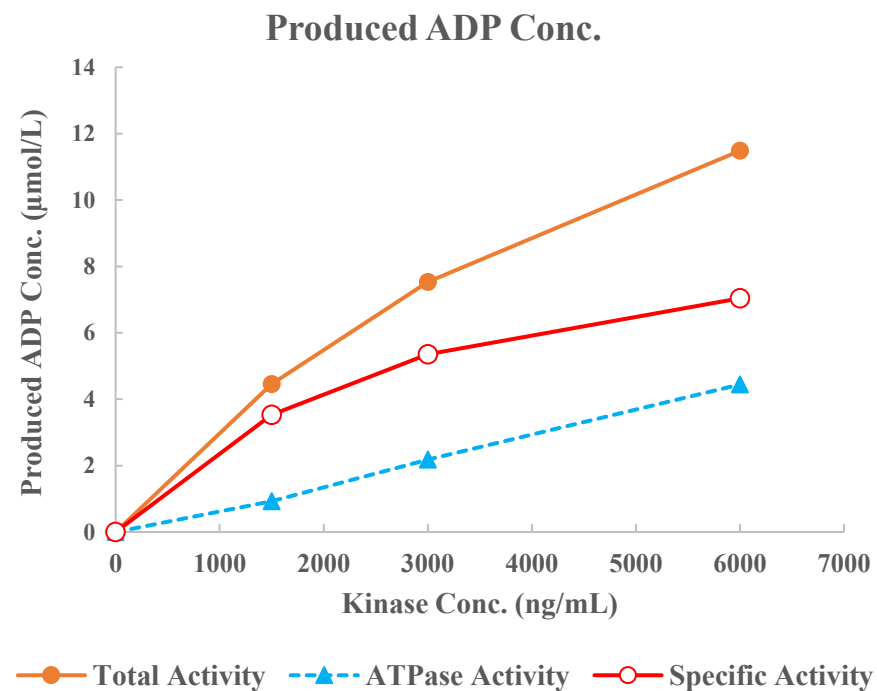
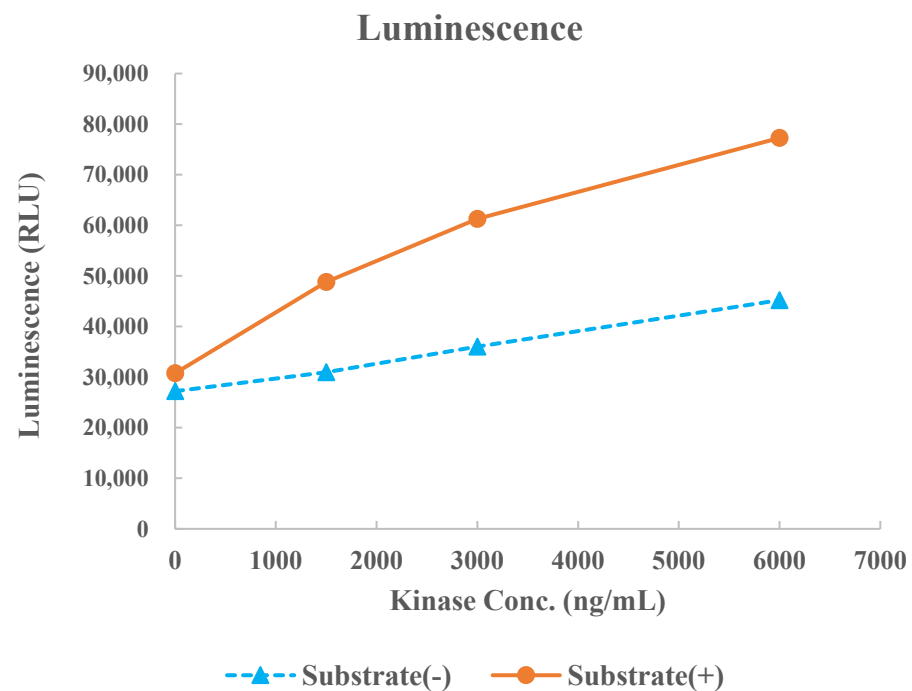


CDK4/CycD1



Kinase	CDK4/CycD1, 04-120
Substrate	Biotin-DYRKtide, BTSU06
Substrate Conc.	10 μmol/L
ATP Conc.	1000 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour