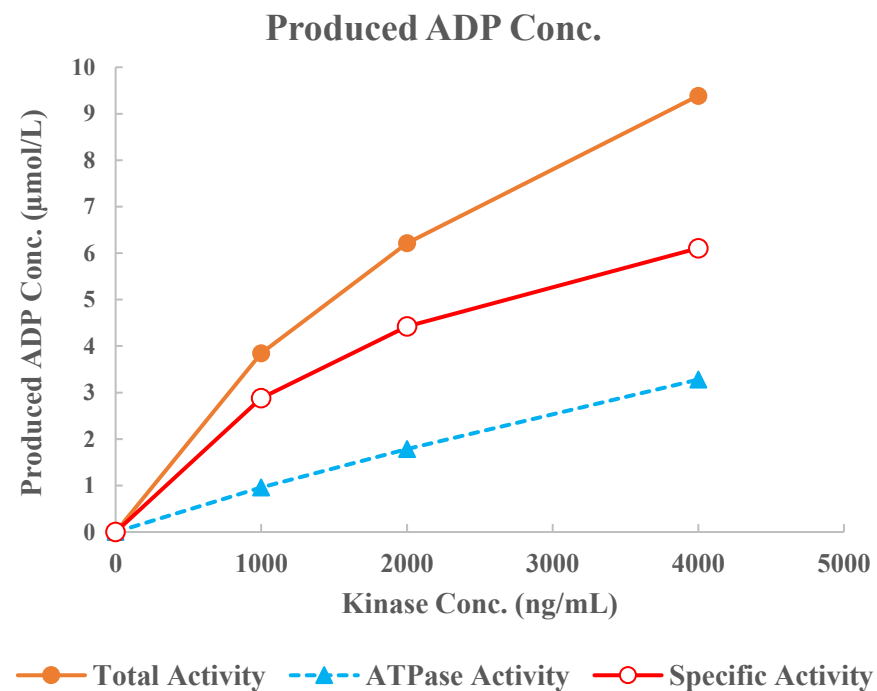
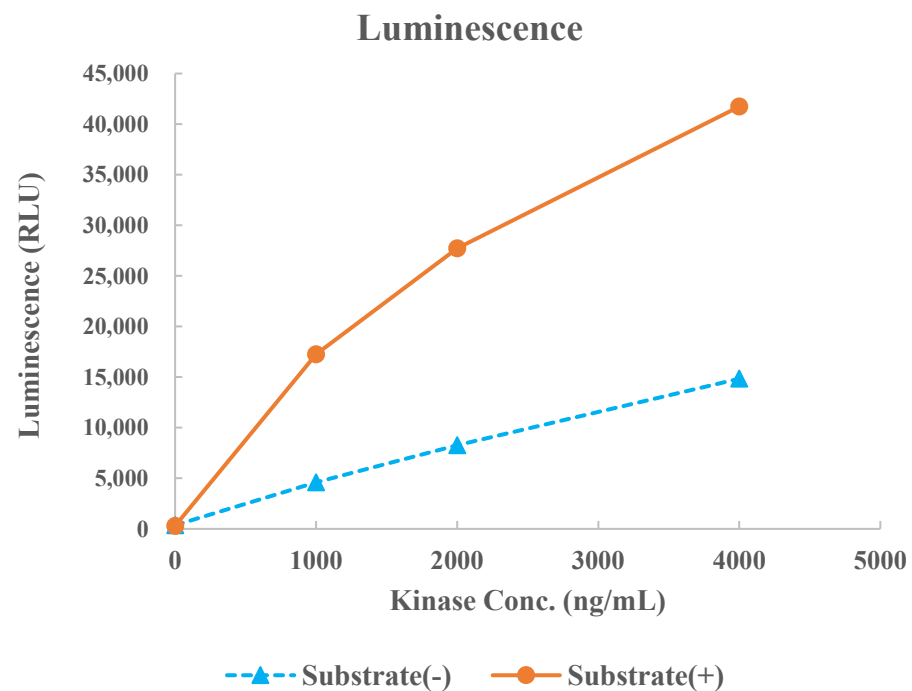


CDK9/CycT1



Kinase	CDK9/CycT1, 04-110
Substrate	Biotin-CDK9 substrate, BTSU29
Substrate Conc.	50 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour