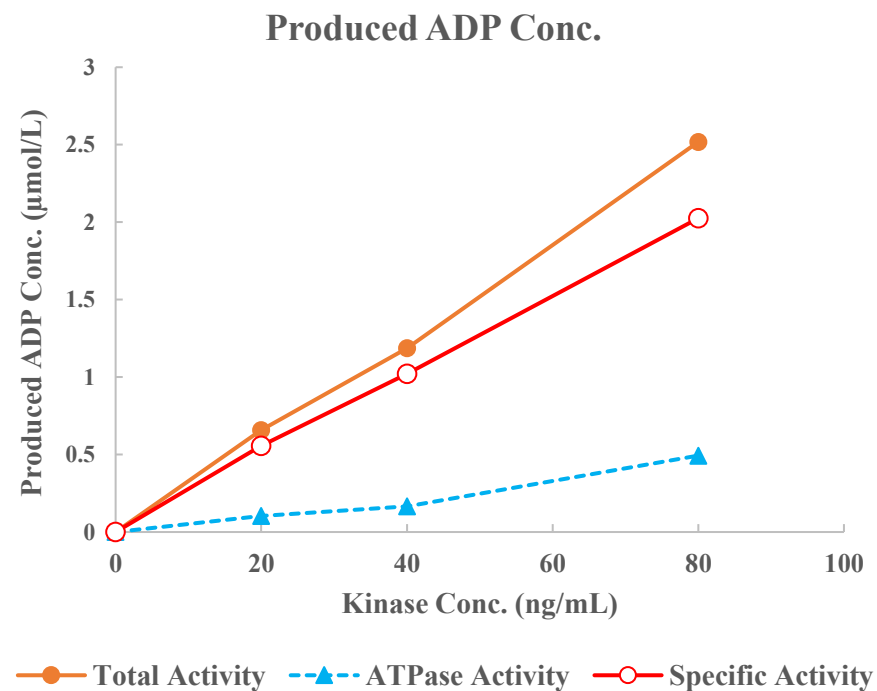
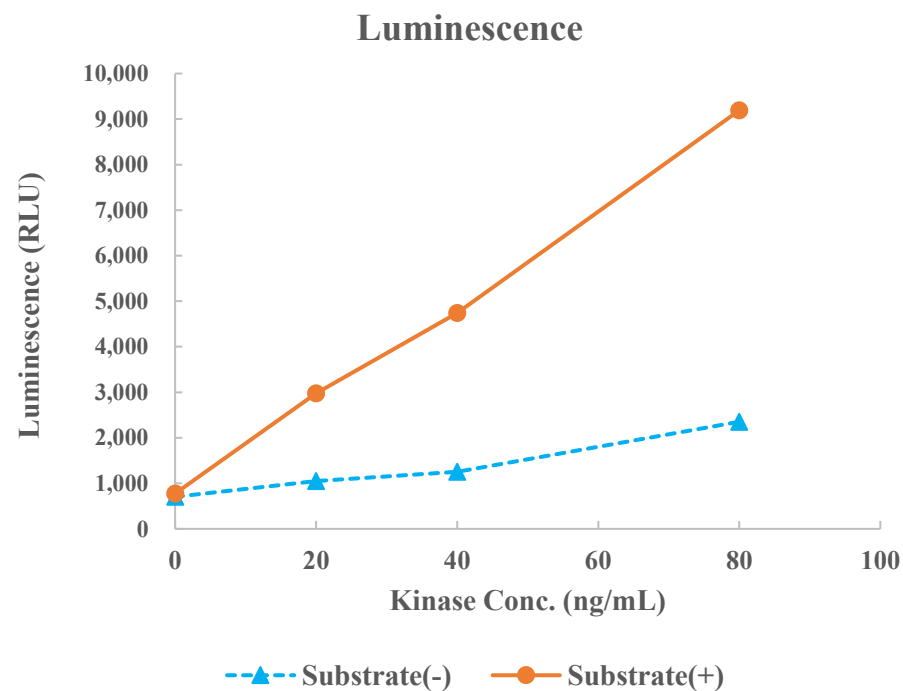


CDK2/CycA2



Kinase	CDK2/CycA2, 04-103
Substrate	Biotin-Modified Histone H1, BTSU13
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour