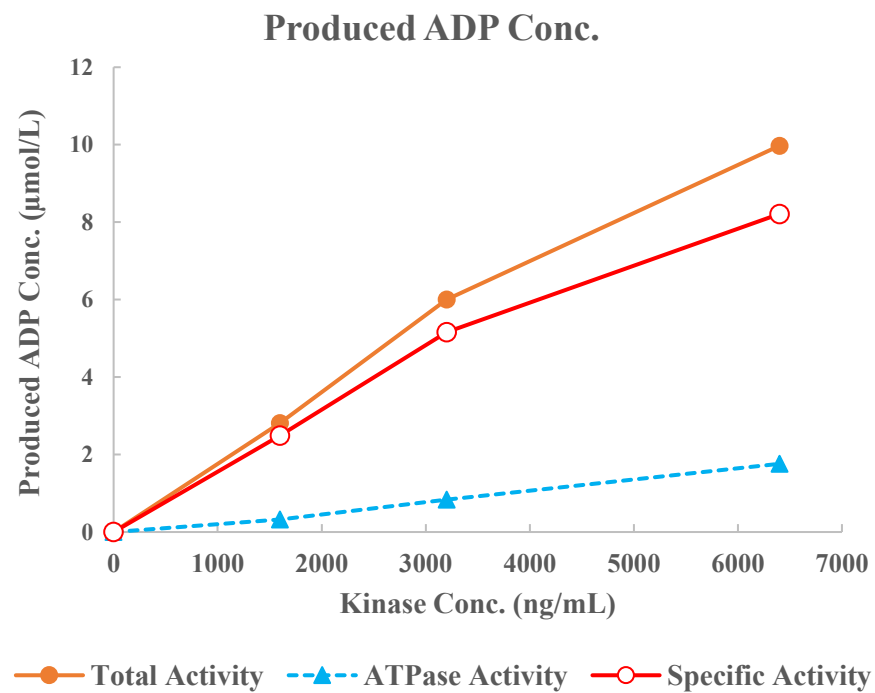
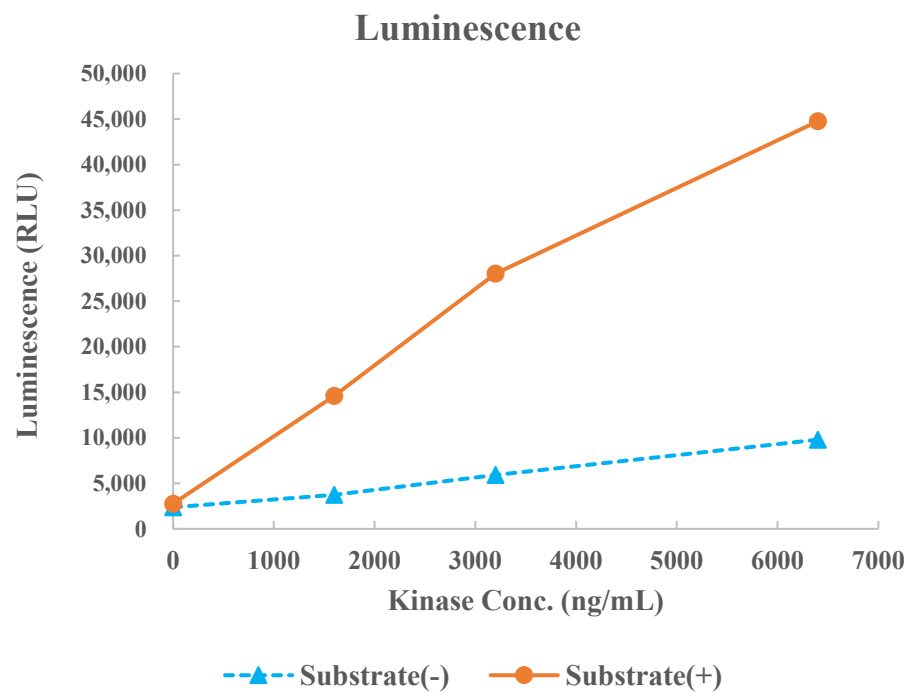


CDK7/CycH/MAT1



Kinase	CDK7/CycH/MAT1, 04-108
Substrate	Biotin-CTD3 peptide, BTSU30
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour