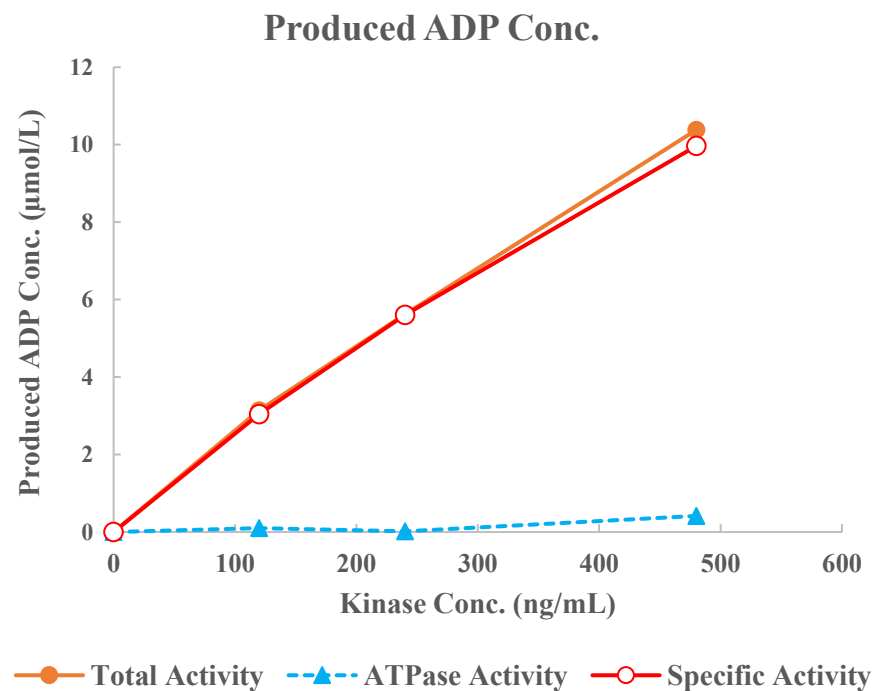
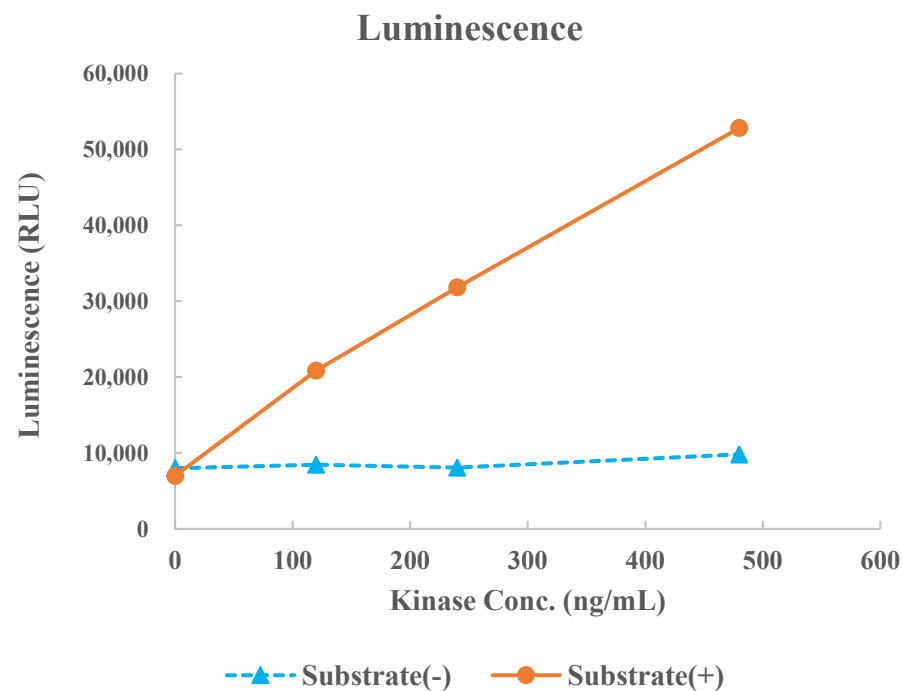


CDK6/CycD3



Kinase	CDK6/CycD3, 04-107
Substrate	Biotin-DYRKtide, BTSU06
Substrate Conc.	50 μmol/L
ATP Conc.	300 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour