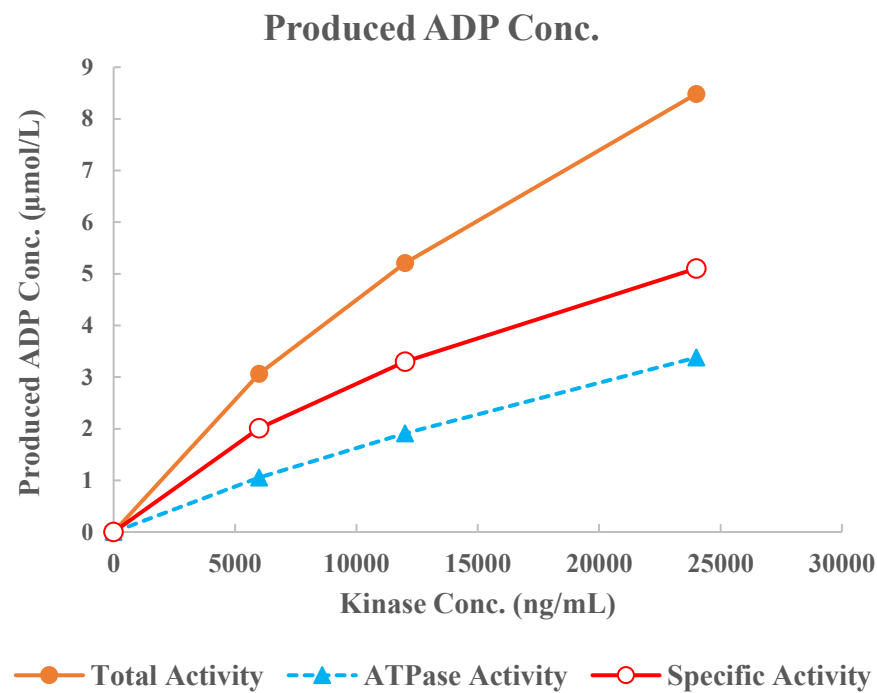
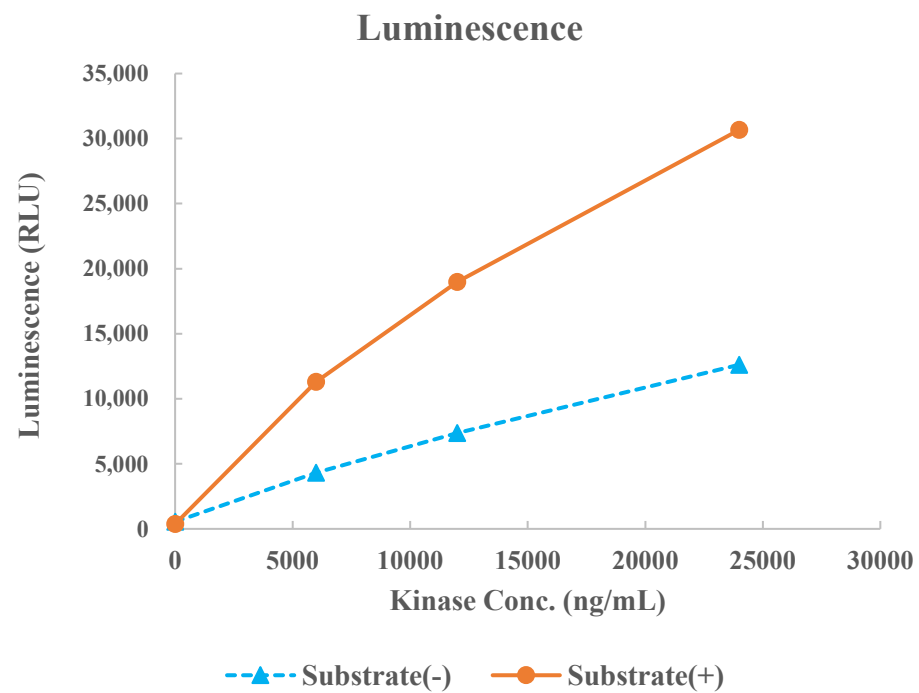


PLK3



Kinase	PLK3, 05-159
Substrate	Biotin-CDC25ctide, BTSU21
Substrate Conc.	50 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour