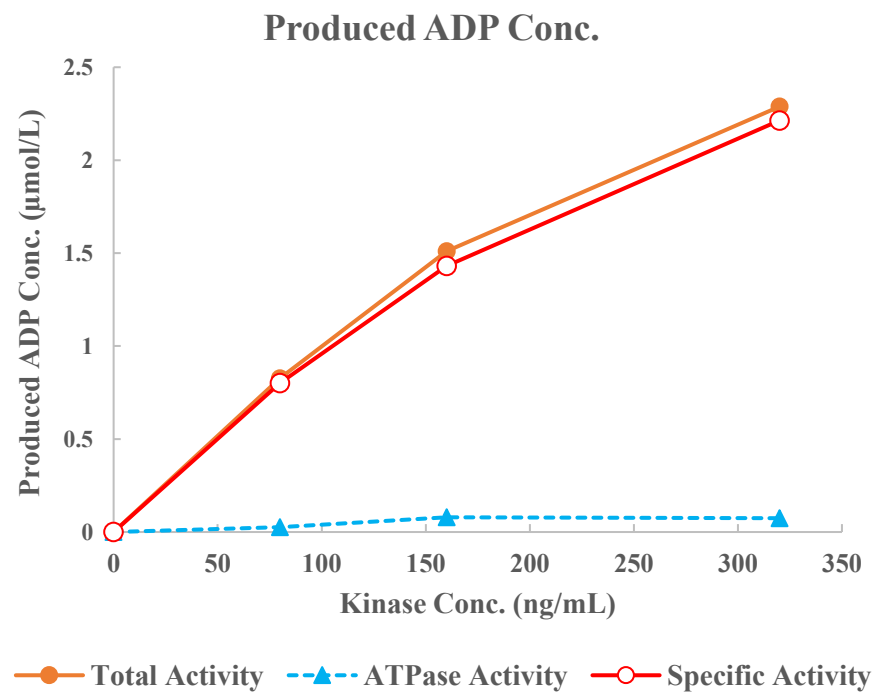
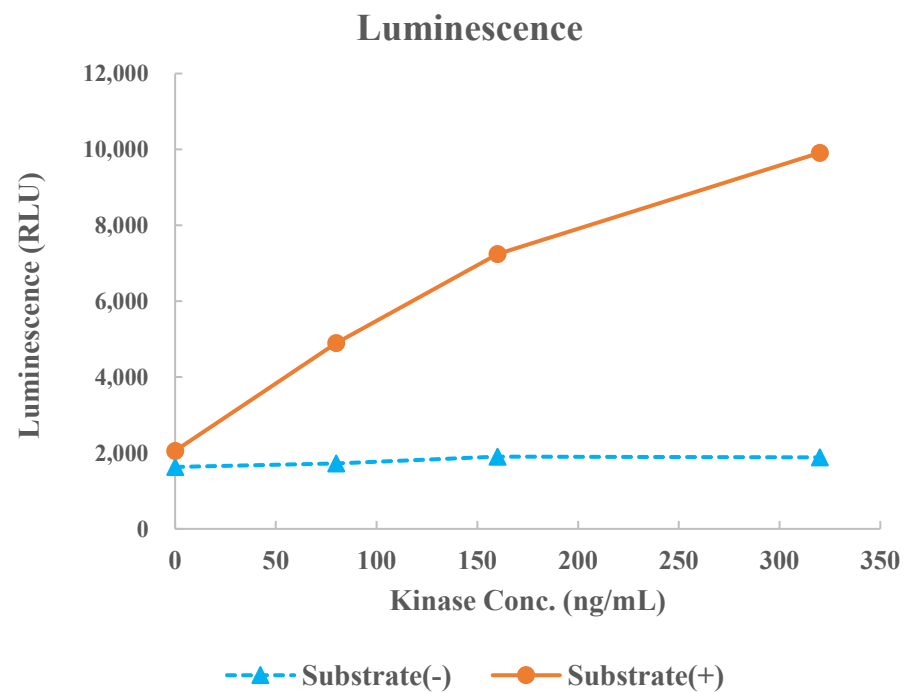


CHK1(CHEK1)



Kinase	CHK1(CHEK1), 02-117
Substrate	Biotin-CHKtide, BTSU08
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour