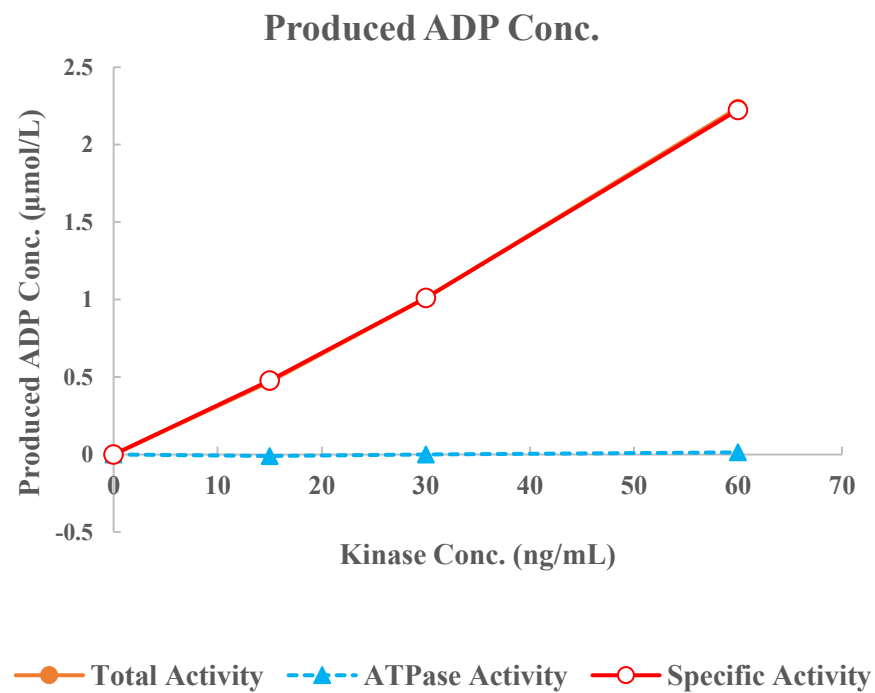
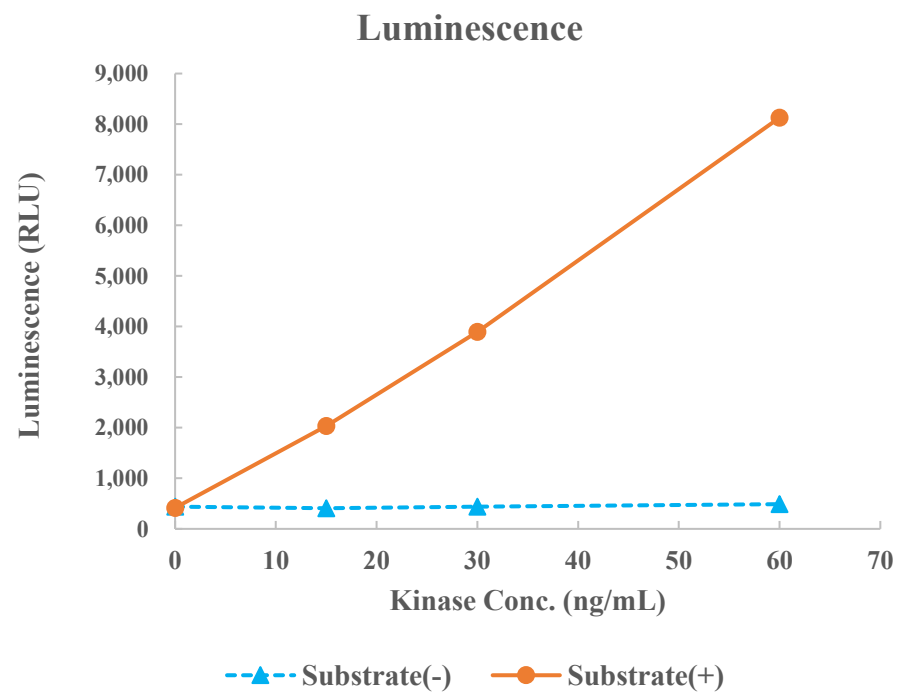


MARK1



Kinase	MARK1, 02-120
Substrate	Biotin-CHKtide, BTSU08
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour