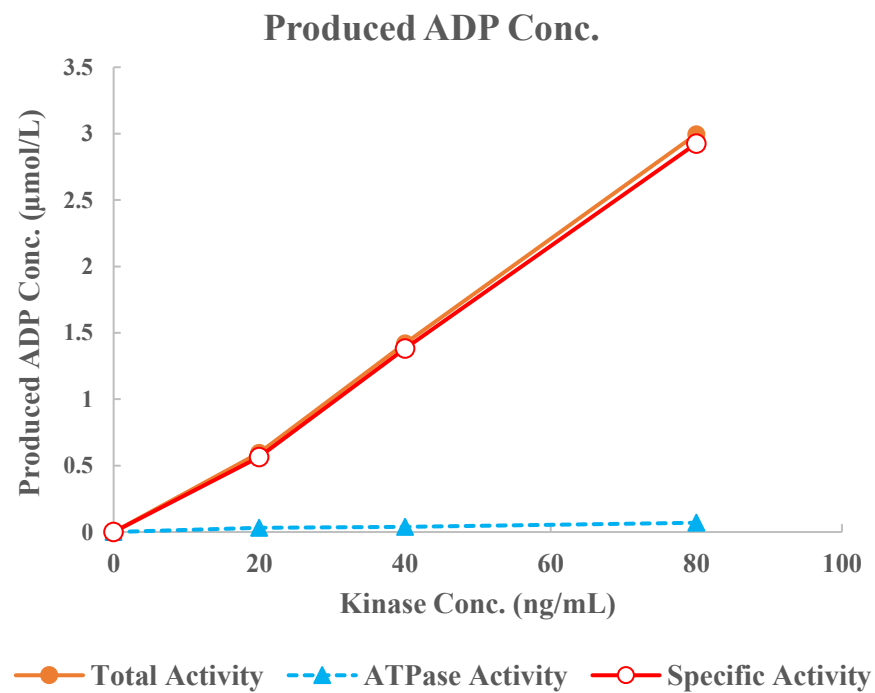
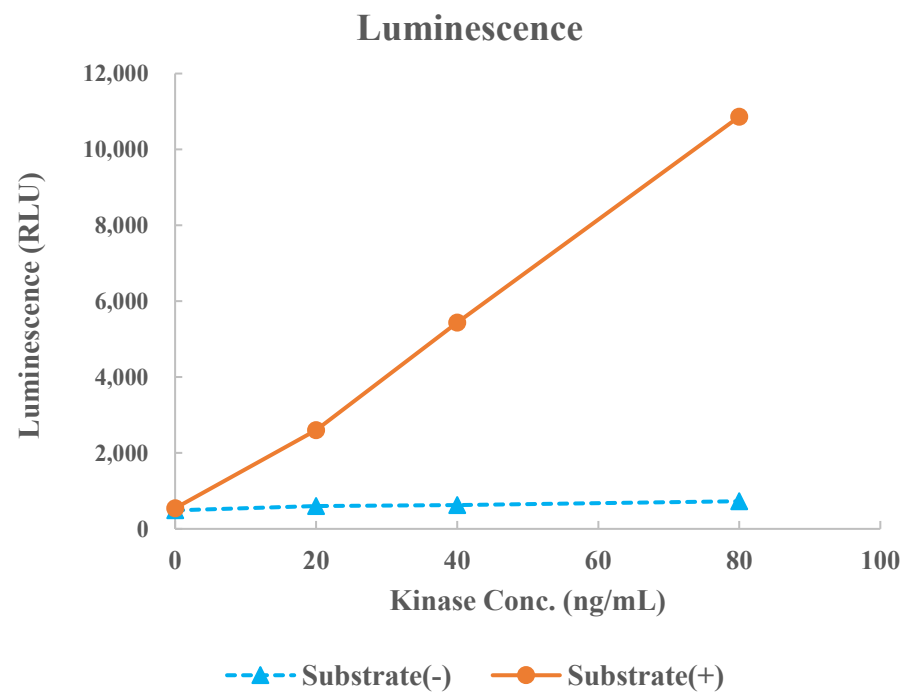


MARK4



Kinase	MARK4, 02-123
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