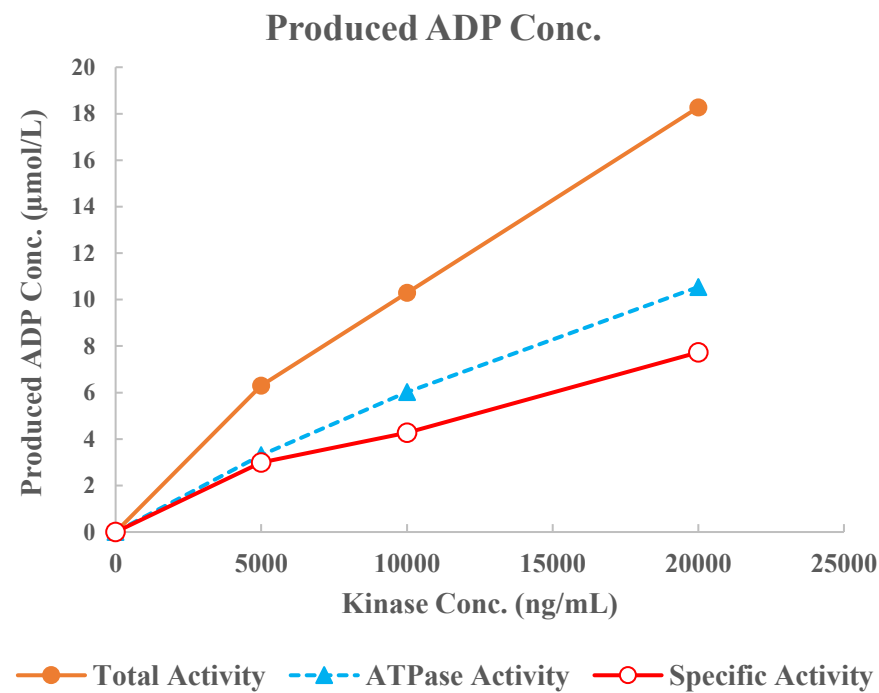
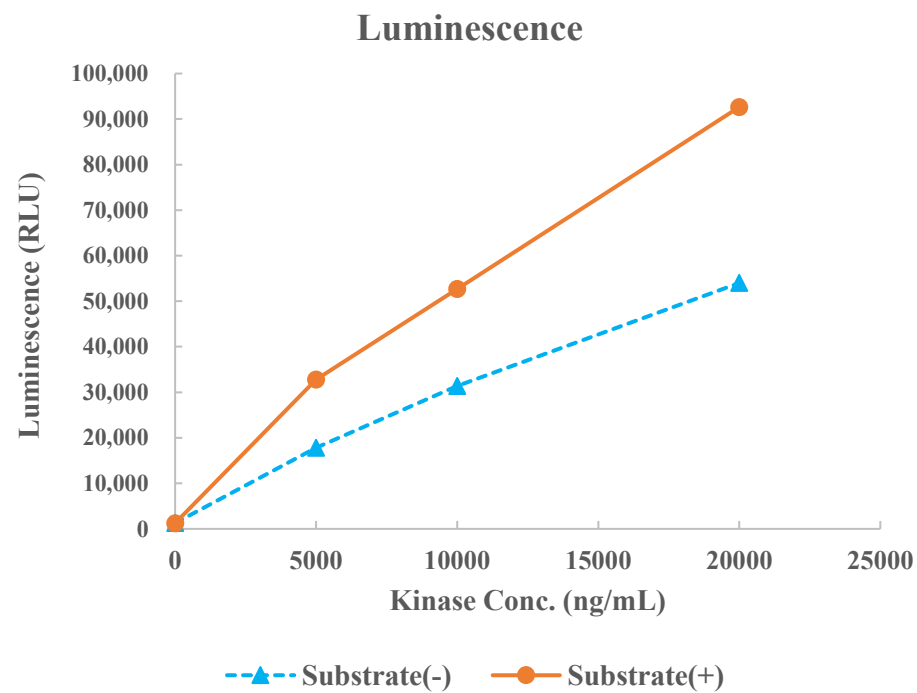


TAOK2



Kinase	TAOK2, 07-133
Substrate	Biotin-TAOKtide, BTSU46
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	10 μmol/L Cantharidin
Reaction Time	1 hour