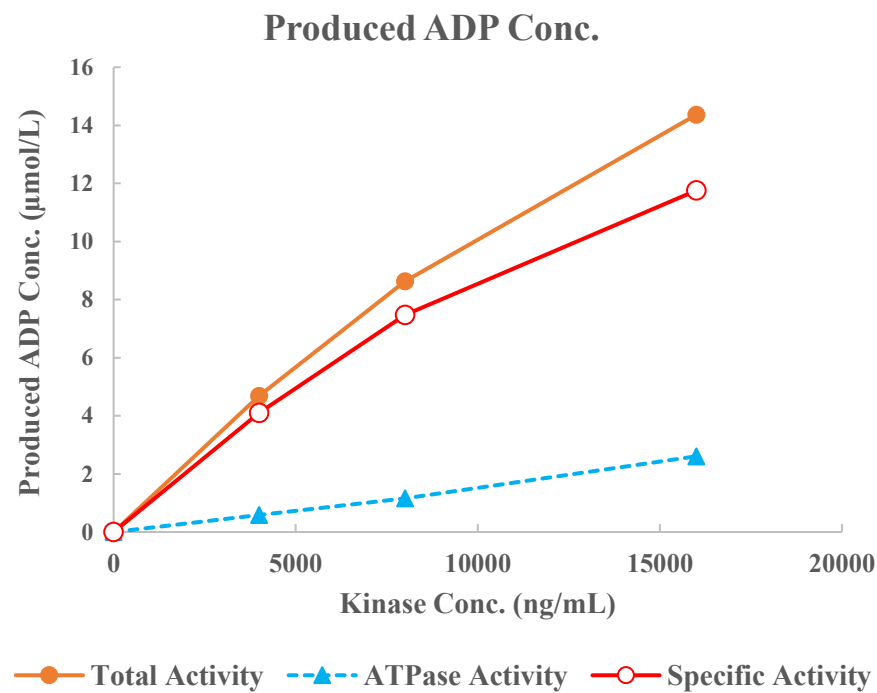
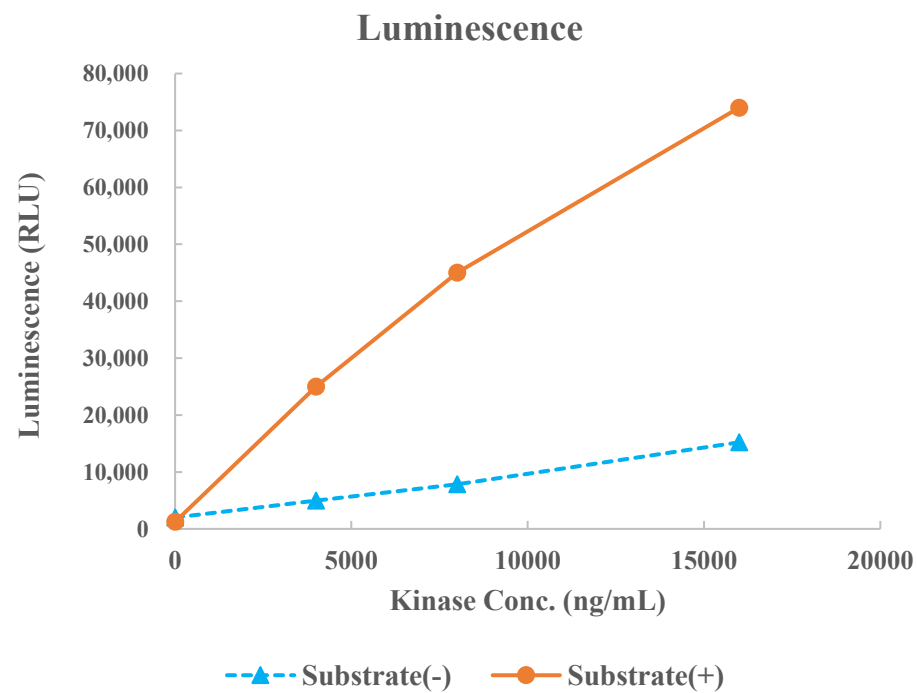


PDHK4(PDK4)



Kinase	PDHK4(PDK4), 10-125
Substrate	Biotin-PDHKtide, BTSU35
Substrate Conc.	50 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg, K
Metal Conc.	5, 25 mmol/L
Additives	-
Reaction Time	1 hour