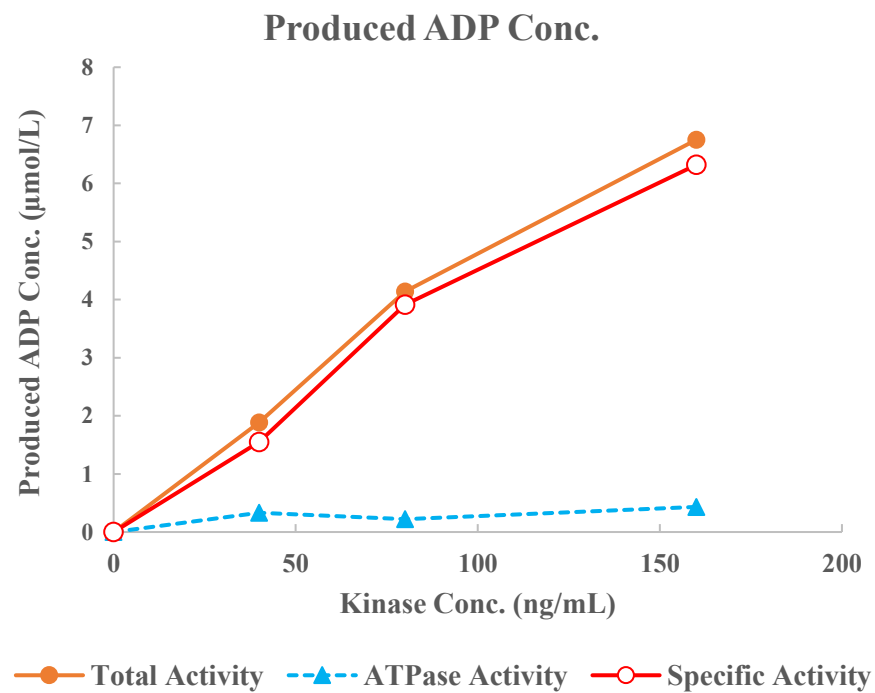
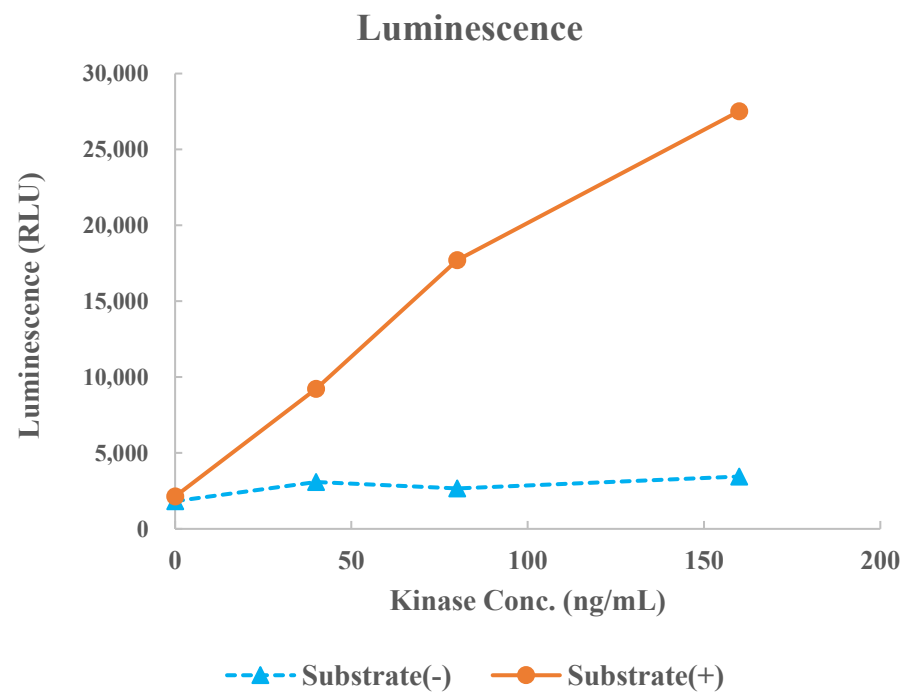


DYRK1B



Kinase	DYRK1B, 04-131
Substrate	Biotin-DYRKtide, BTSU06
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour