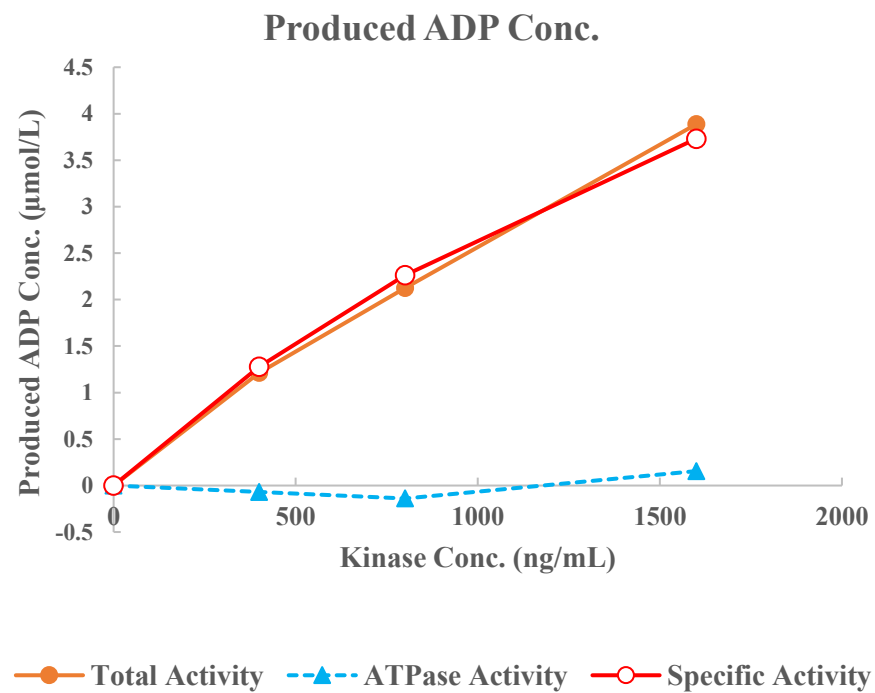
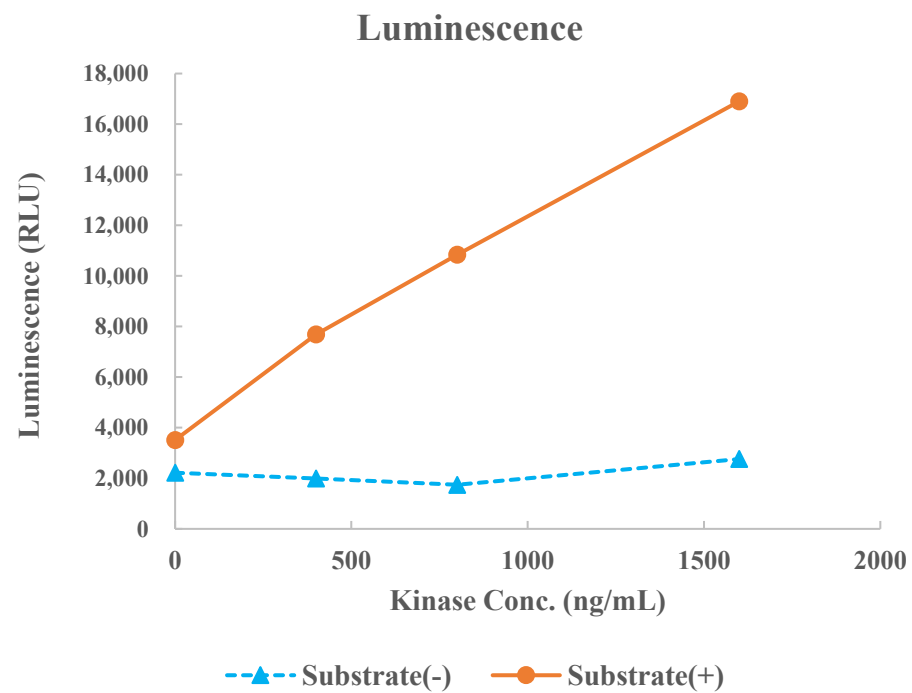


BRSK2



Kinase	BRSK2, 02-116
Substrate	Biotin-CHKtide, BTSU08
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour