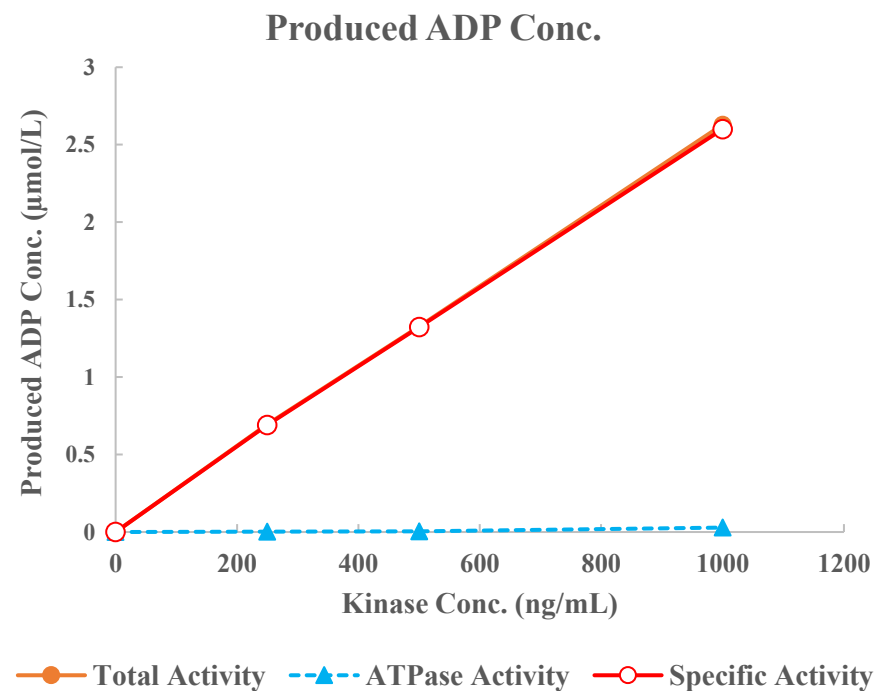
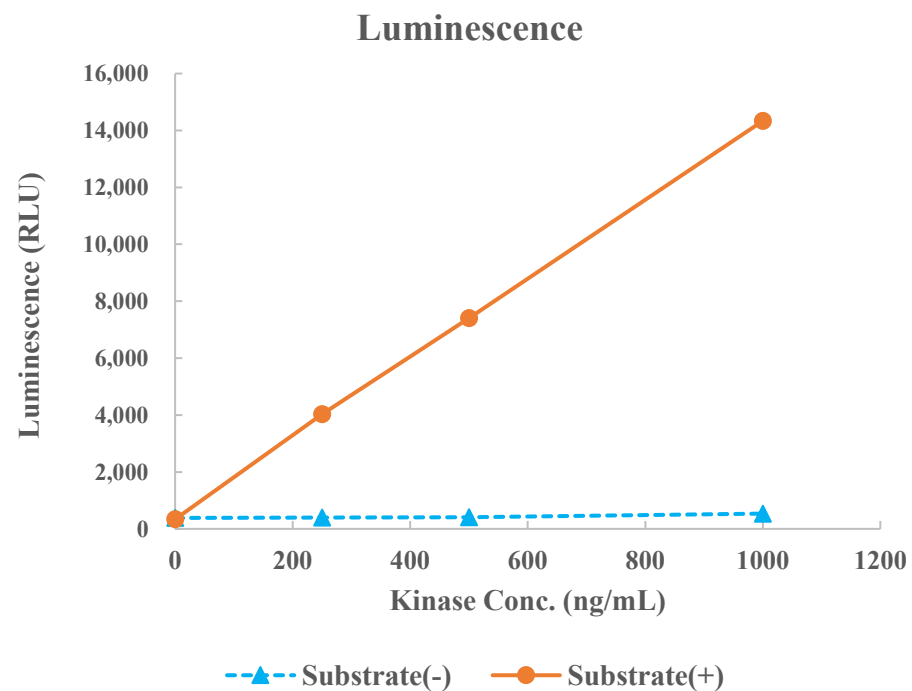


TSSK2



Kinase	TSSK2, 02-165
Substrate	Biotin-GS peptide, BTSU05
Substrate Conc.	50 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour