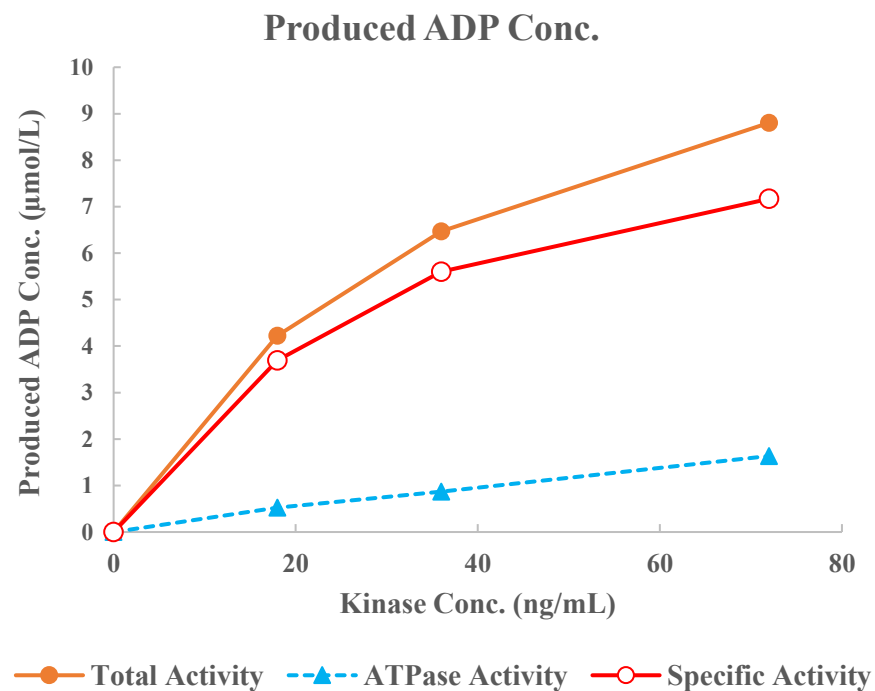
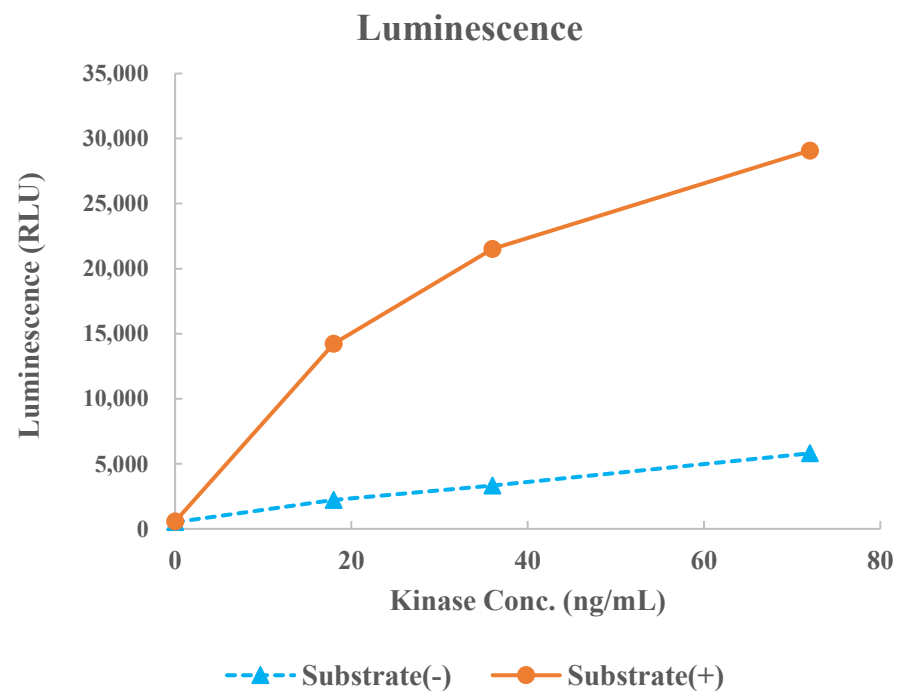


JAK2



Kinase	JAK2, 08-045
Substrate	Biotin-Src tide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour