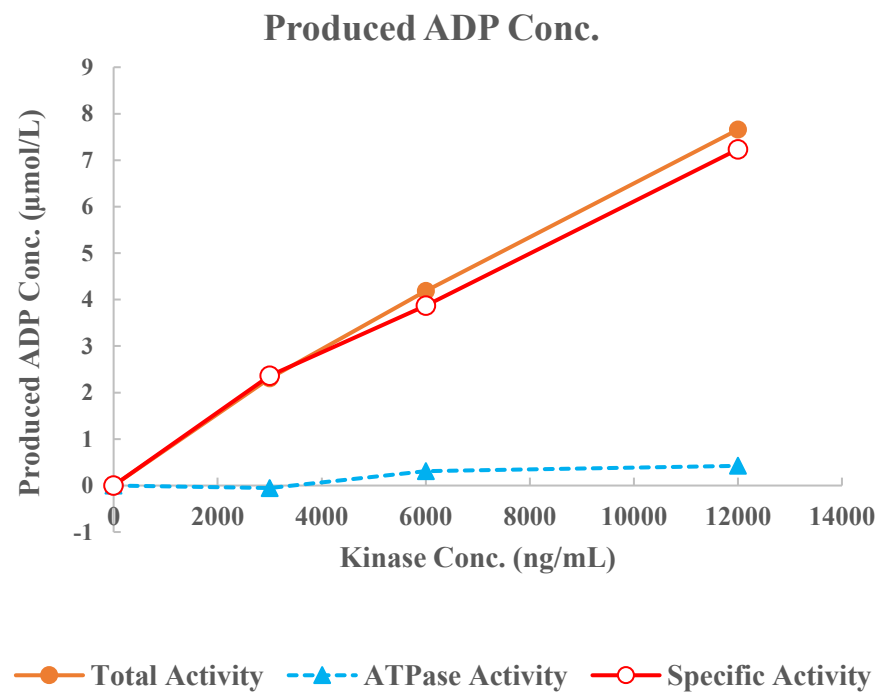
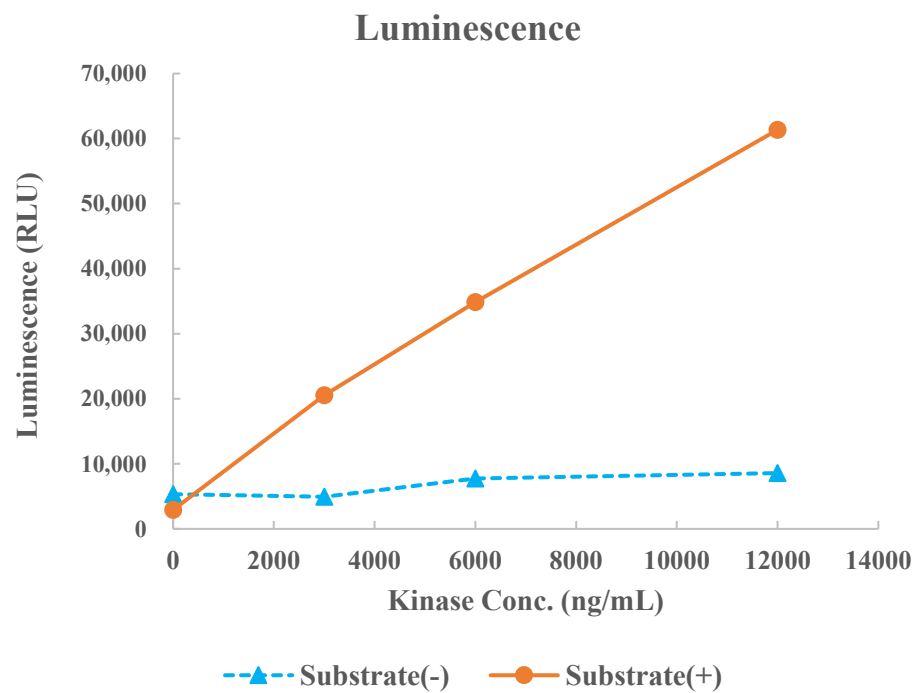


JAK1



Kinase	JAK1, 08-144
Substrate	Biotin-JAK1 substrate peptide, BTSU32
Substrate Conc.	50 μmol/L
ATP Conc.	75 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	25 μmol/L Sodium orthovanadate
Reaction Time	1 hour