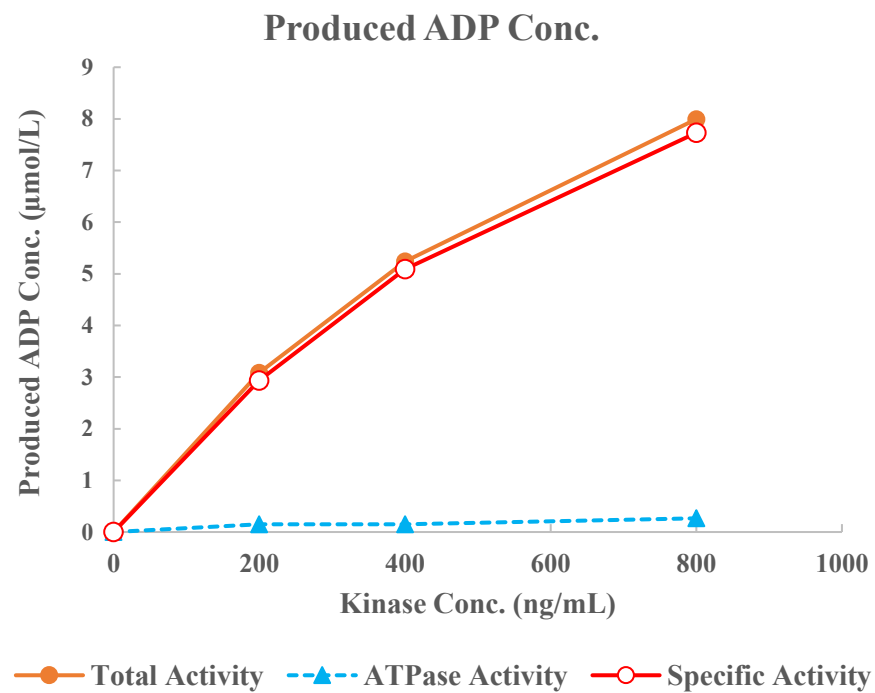
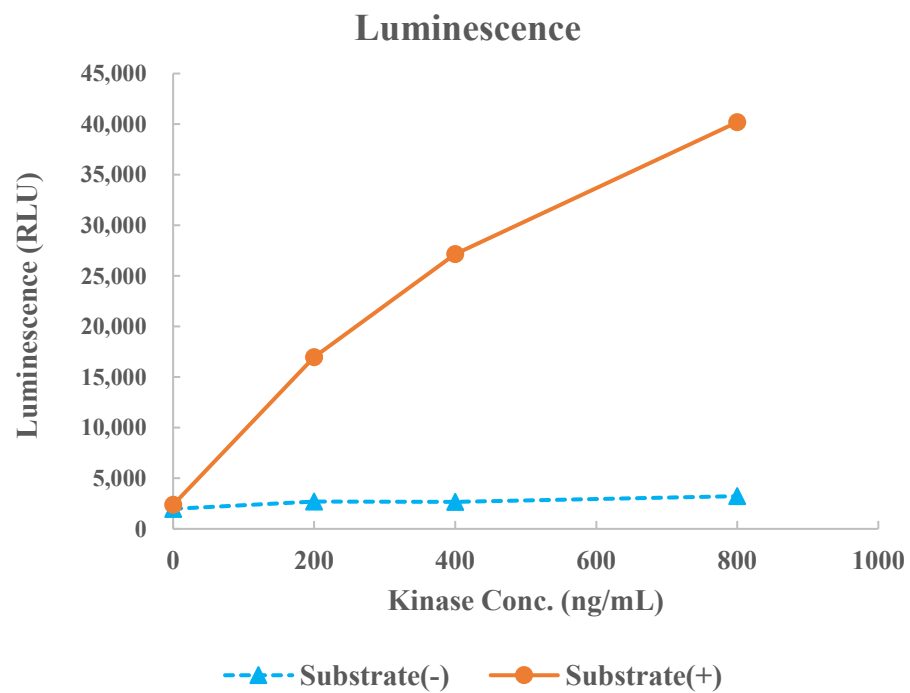


KIT [D816E]



Kinase	KIT [D816E], 08-541
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	25 μmol/L Sodium orthovanadate
Reaction Time	1 hour