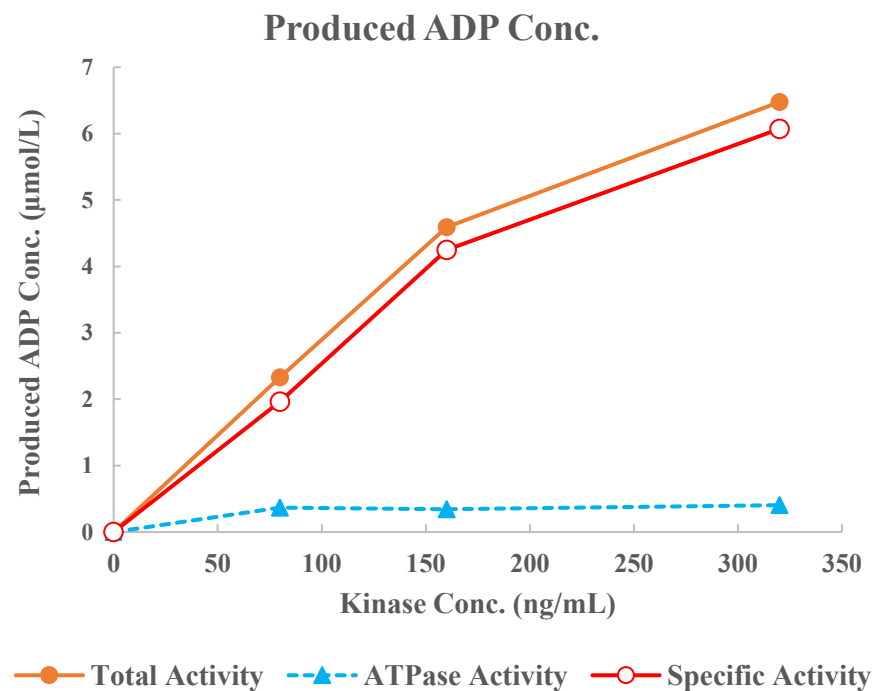
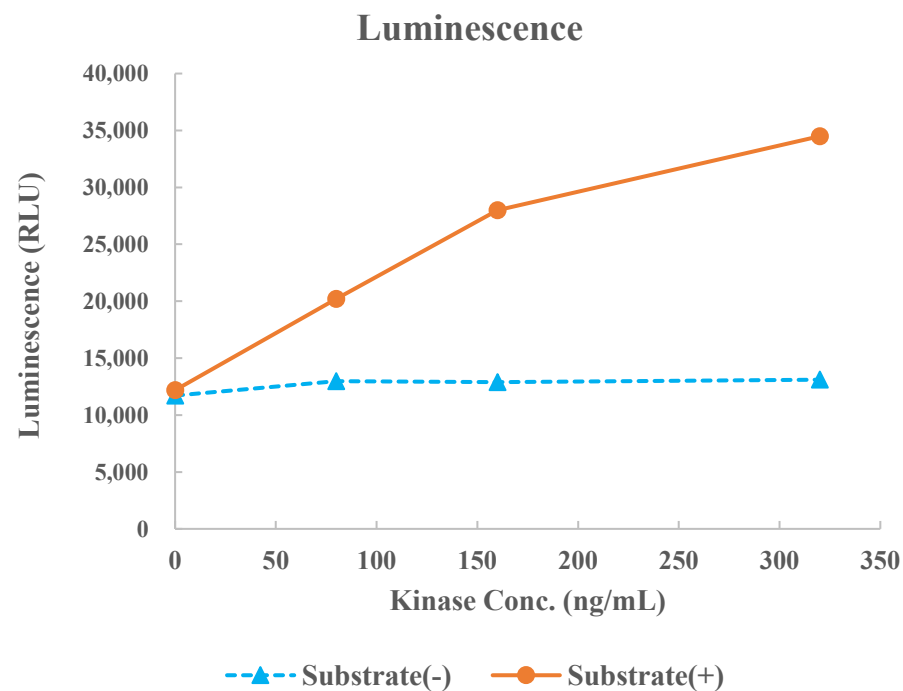


KIT



Kinase	KIT, 08-156
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	400 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	25 μmol/L Sodium orthovanadate
Reaction Time	1 hour