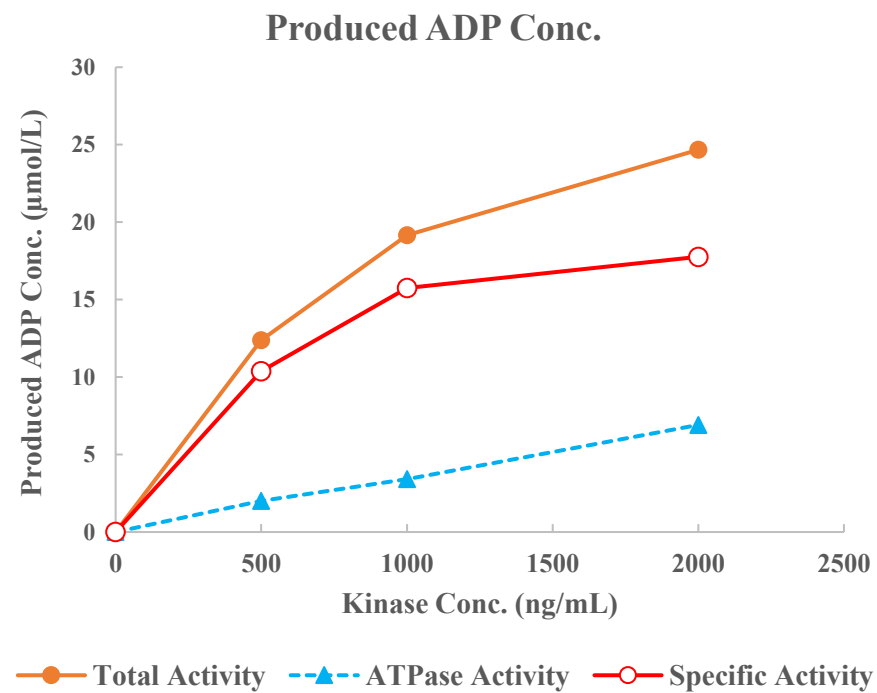
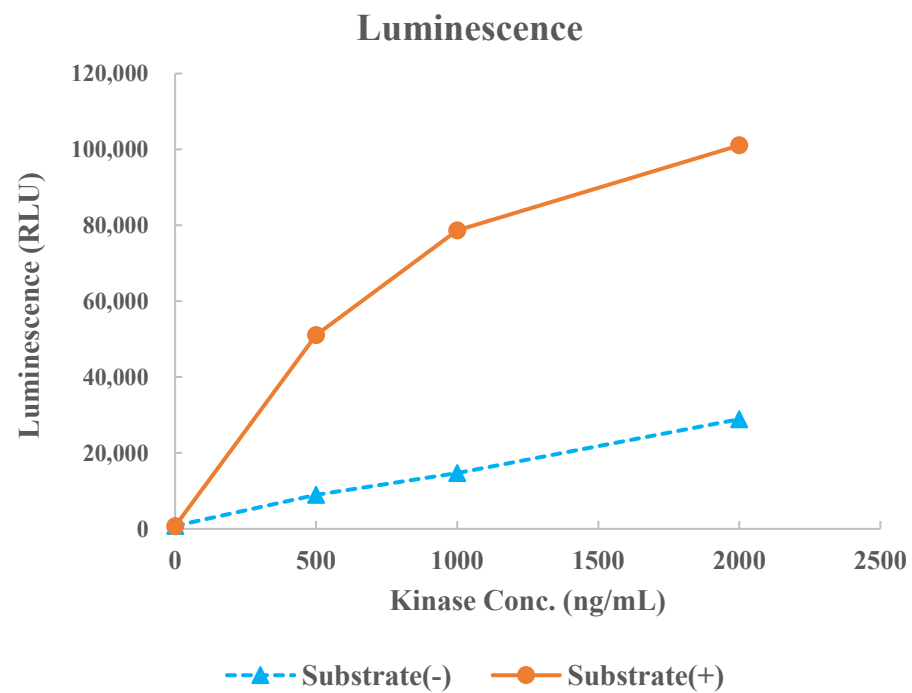


FER



Kinase	FER, 08-139
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	50 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour