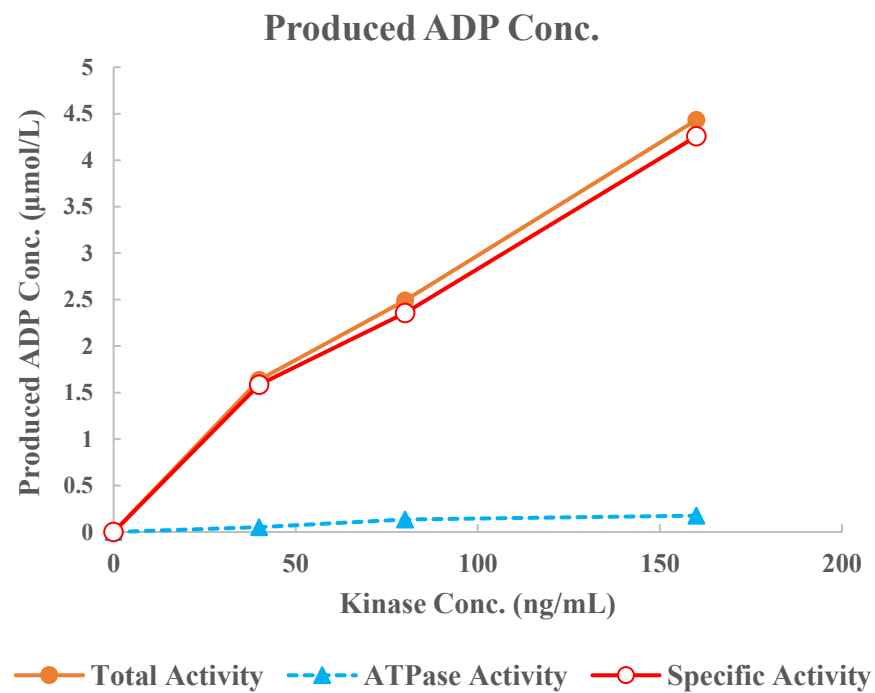
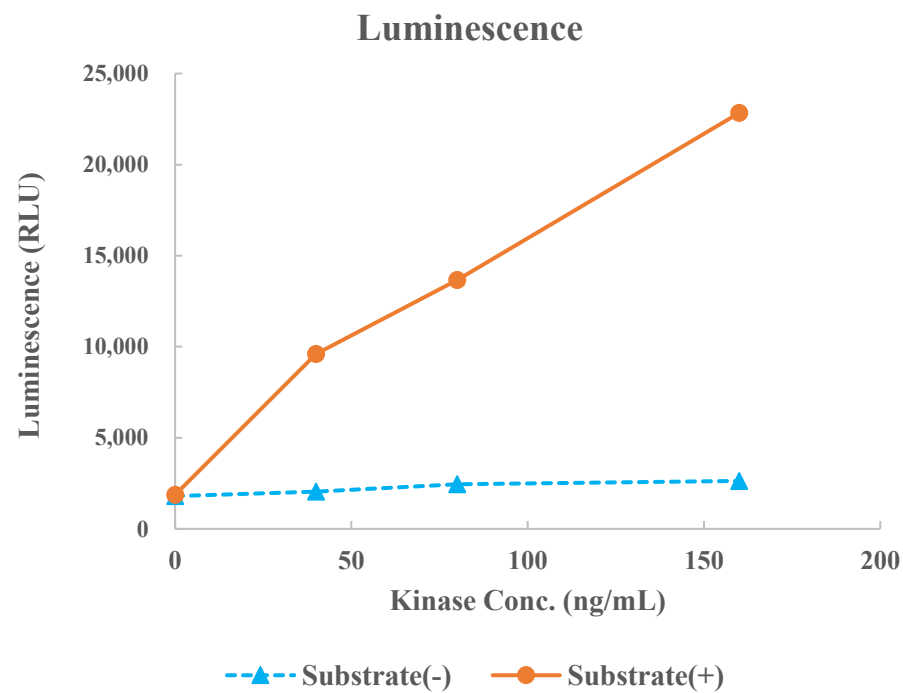


FYN [isoform a]



Kinase	FYN [isoform a], 08-168
Substrate	Biotin-Srcide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour