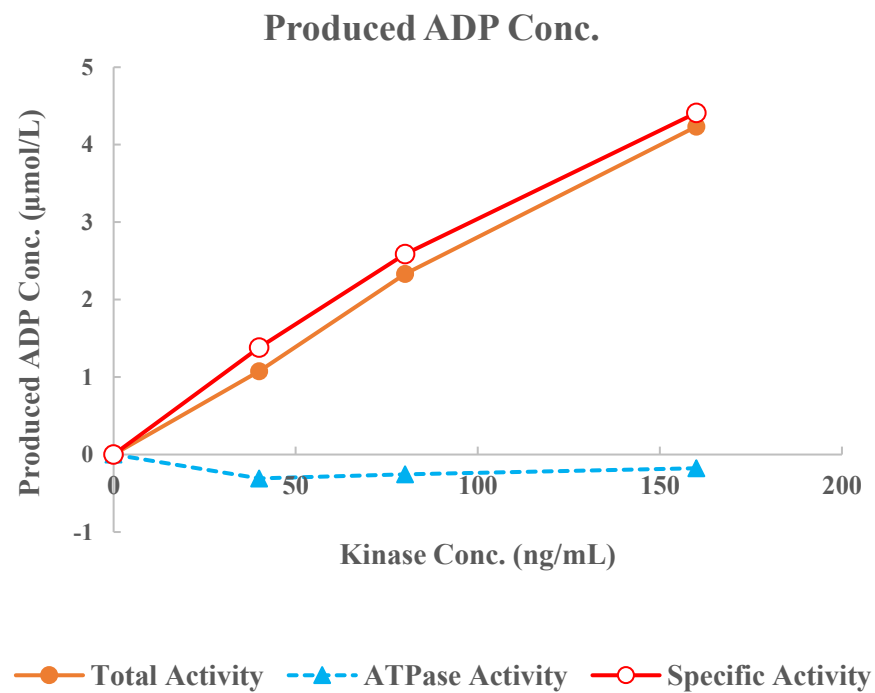
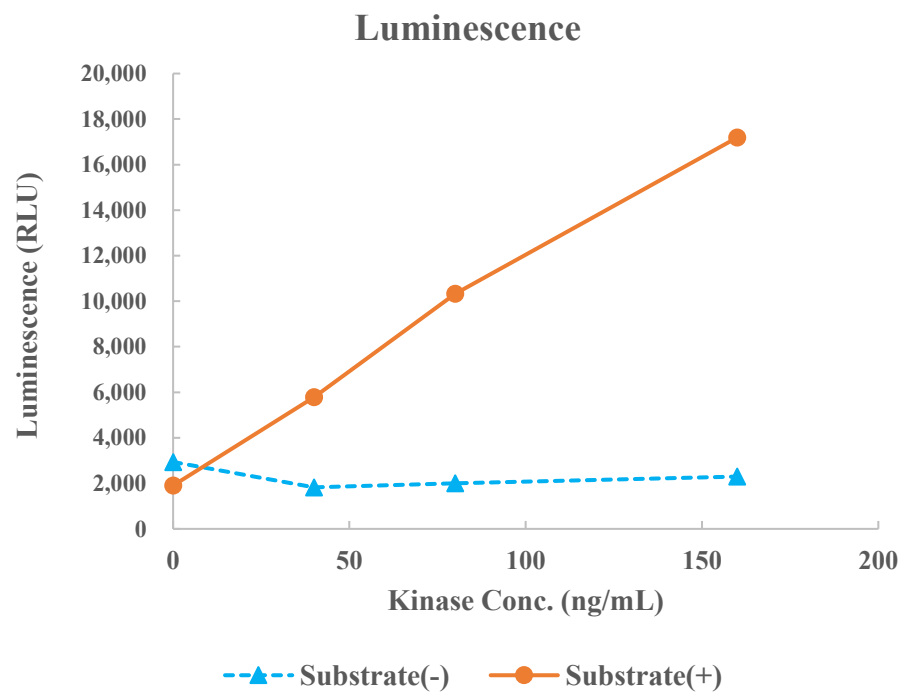


FLT3



Kinase	FLT3, 08-154
Substrate	Biotin-Srcptide, BTSU01
Substrate Conc.	10 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour