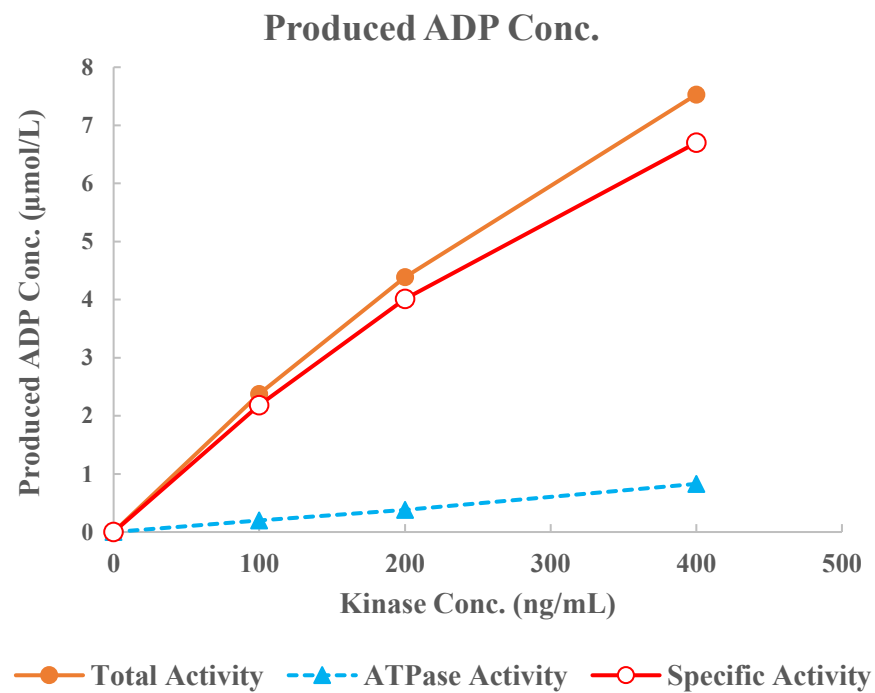
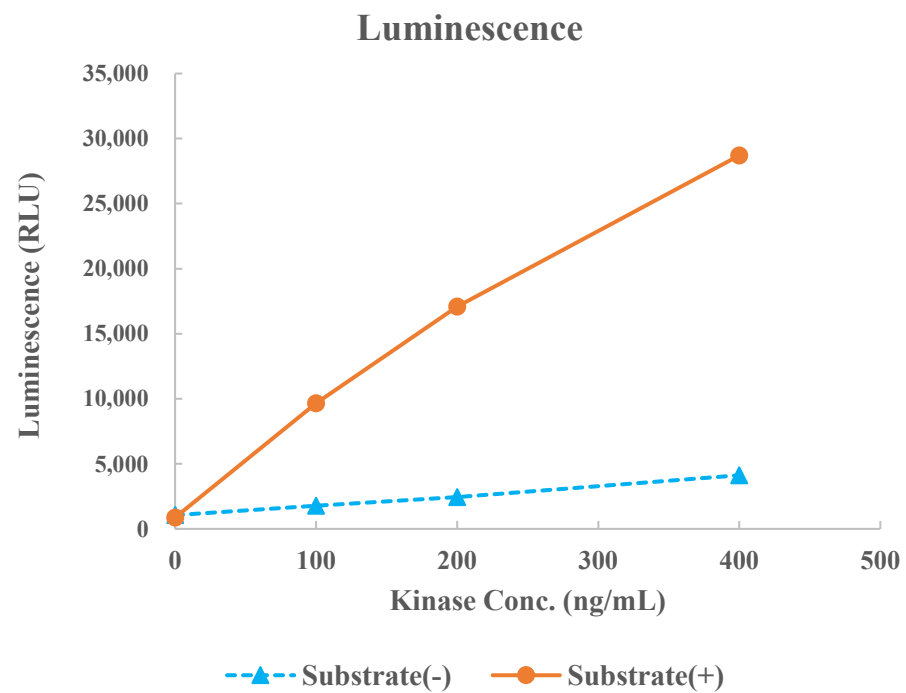


ABL(ABL1)



Kinase	ABL(ABL1), 08-001
Substrate	Biotin-ABLtide, BTSU16
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour