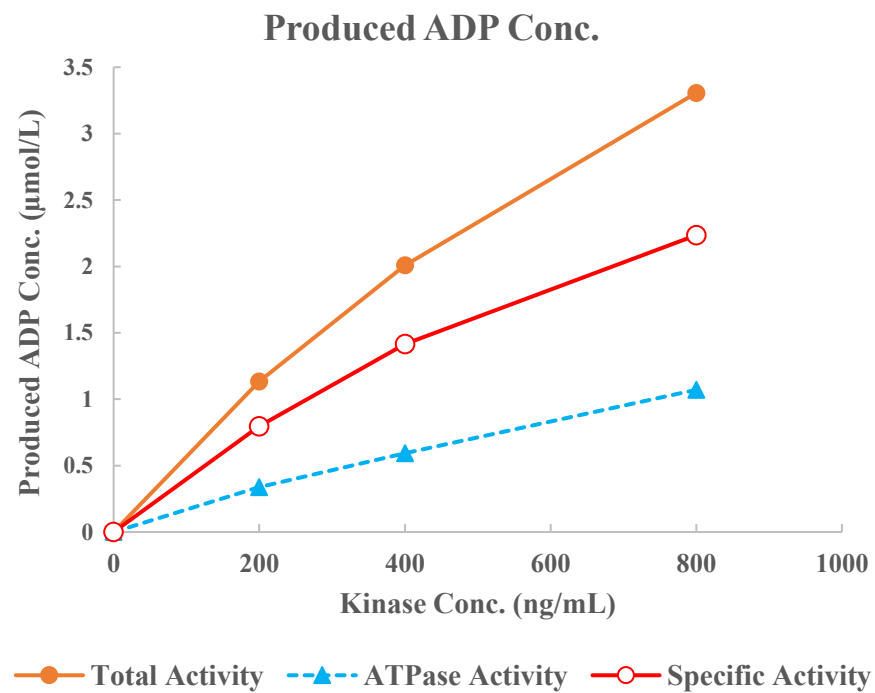
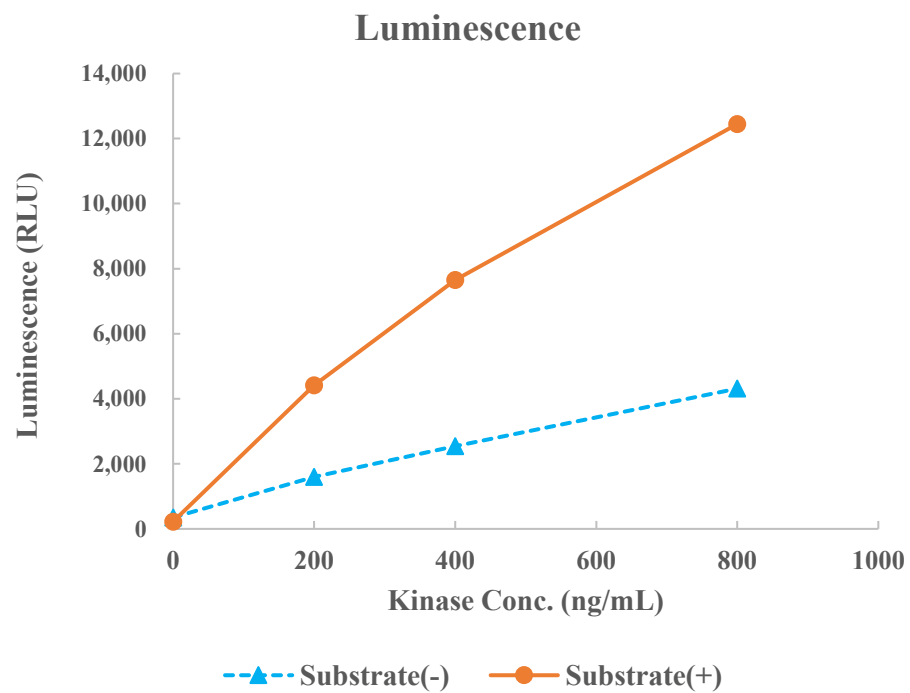


ABL(ABL1) [T315I]



Kinase	ABL(ABL1) [T315I], 08-093
Substrate	Biotin-ABLtide, BTSU16
Substrate Conc.	10 μmol/L
ATP Conc.	5 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour