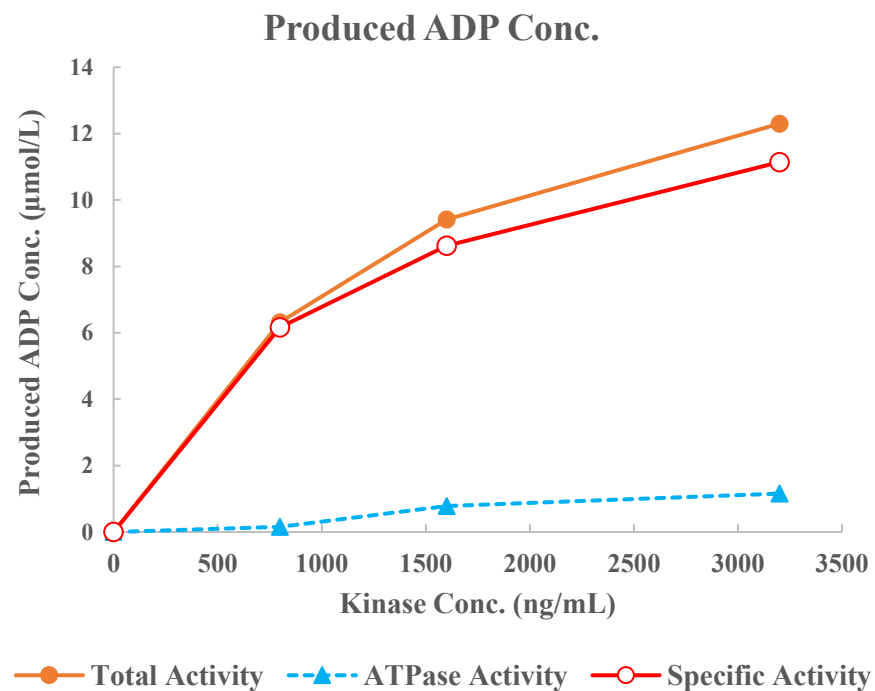
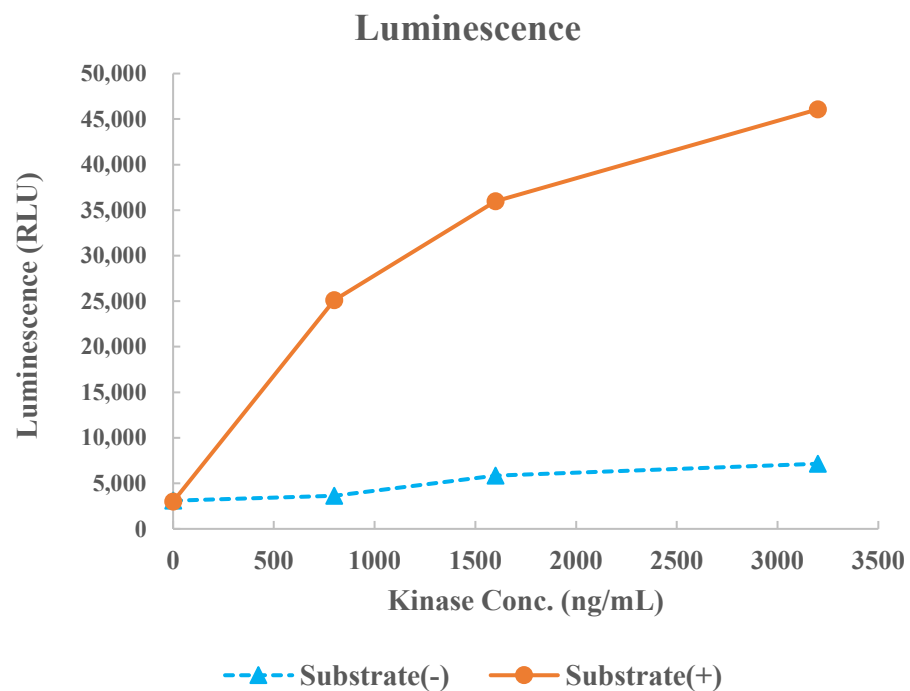


MST3(STK24)



Kinase	MST3(STK24), 07-118
Substrate	Biotin-Moesin-derived peptide, BTSU18
Substrate Conc.	10 μmol/L
ATP Conc.	75 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour