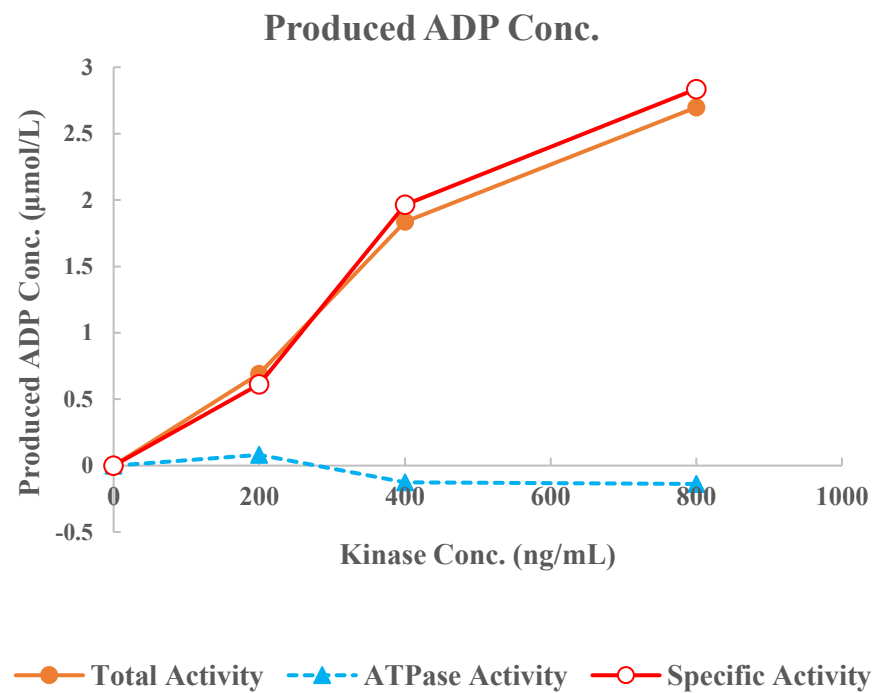
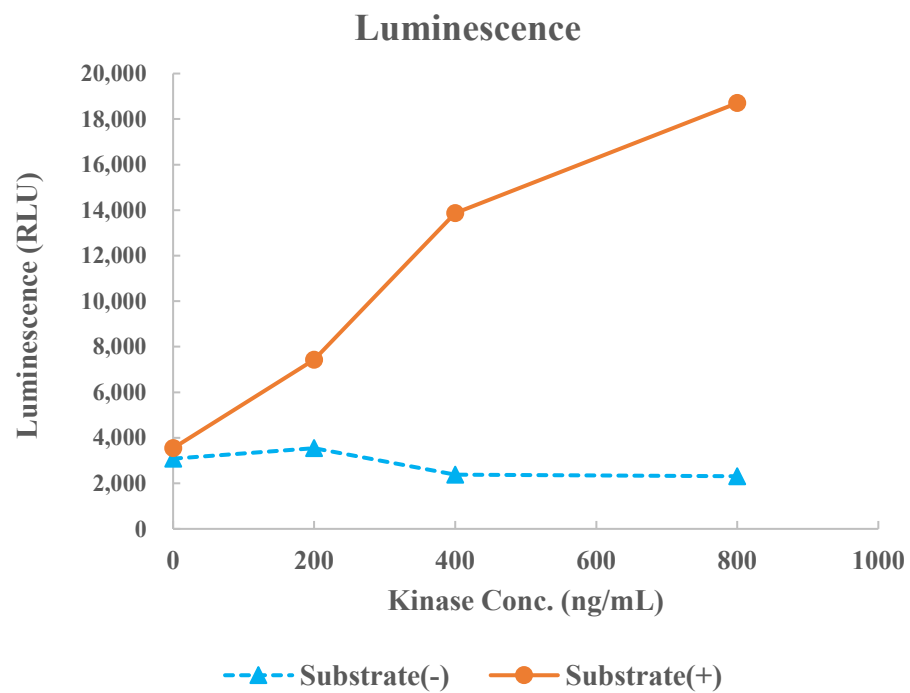


LOK(STK10)



Kinase	LOK(STK10), 07-115
Substrate	Biotin-Moesin-derived peptide, BTSU18
Substrate Conc.	10 μmol/L
ATP Conc.	100 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour