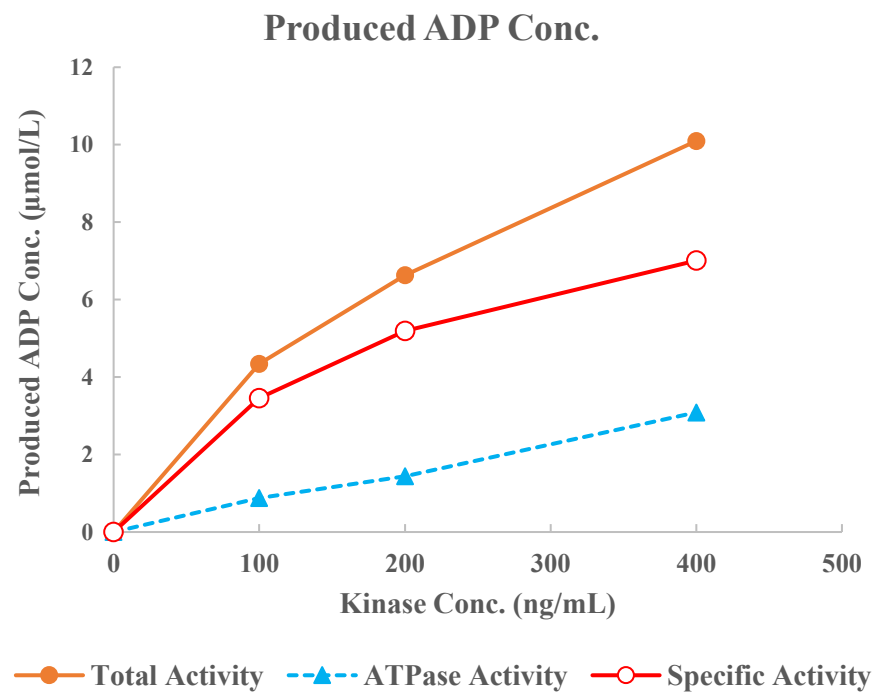
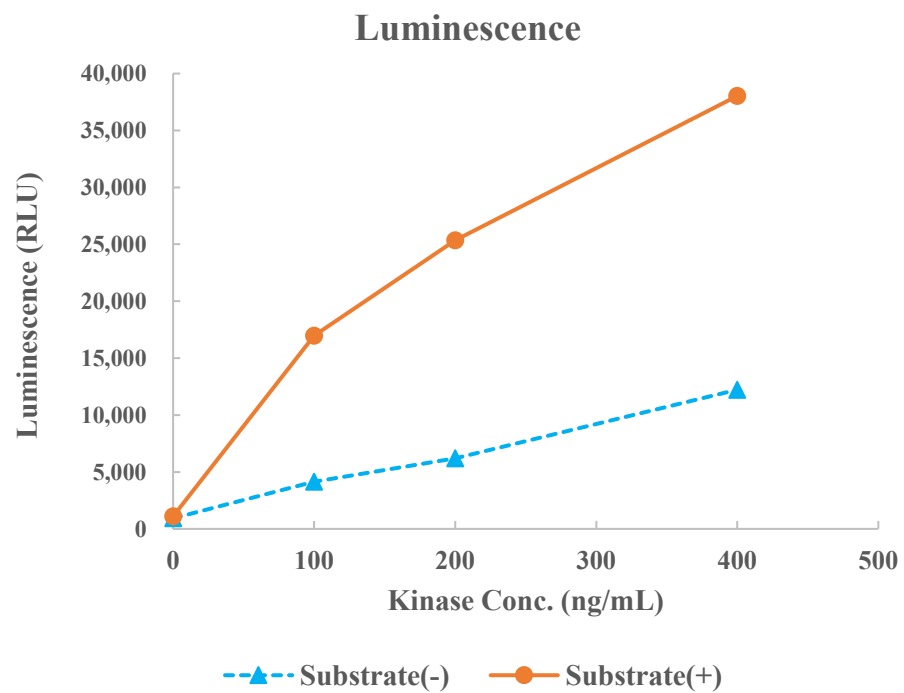


AurA(AURKA)



Kinase	AurA(AURKA), 05-101
Substrate	Biotin-Kemptide, BTSU09
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour