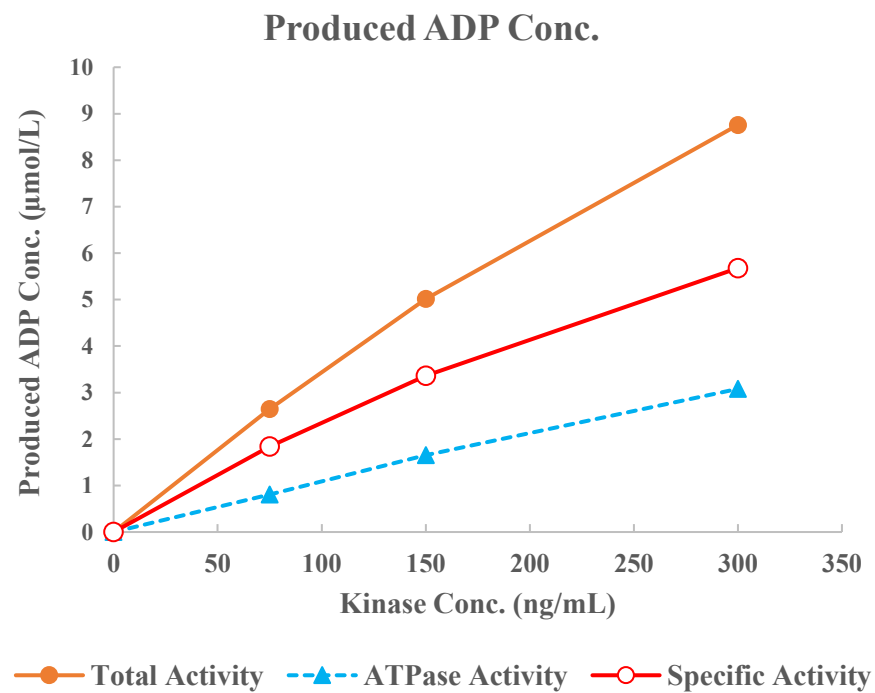
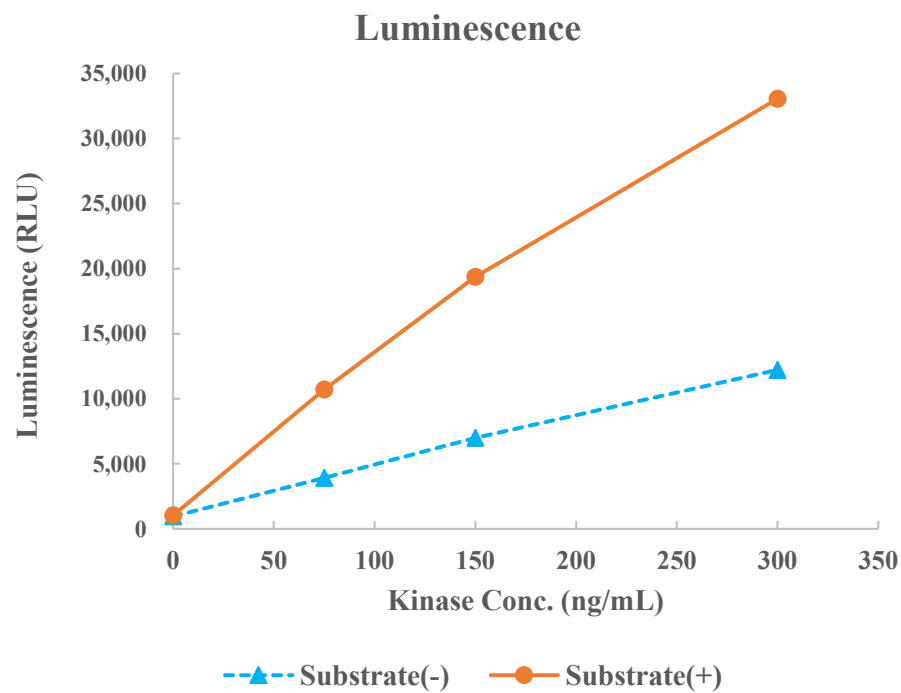


AurB(AURKB)/INCENP



Kinase	AurB(AURKB)/INCENP, 05-102
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