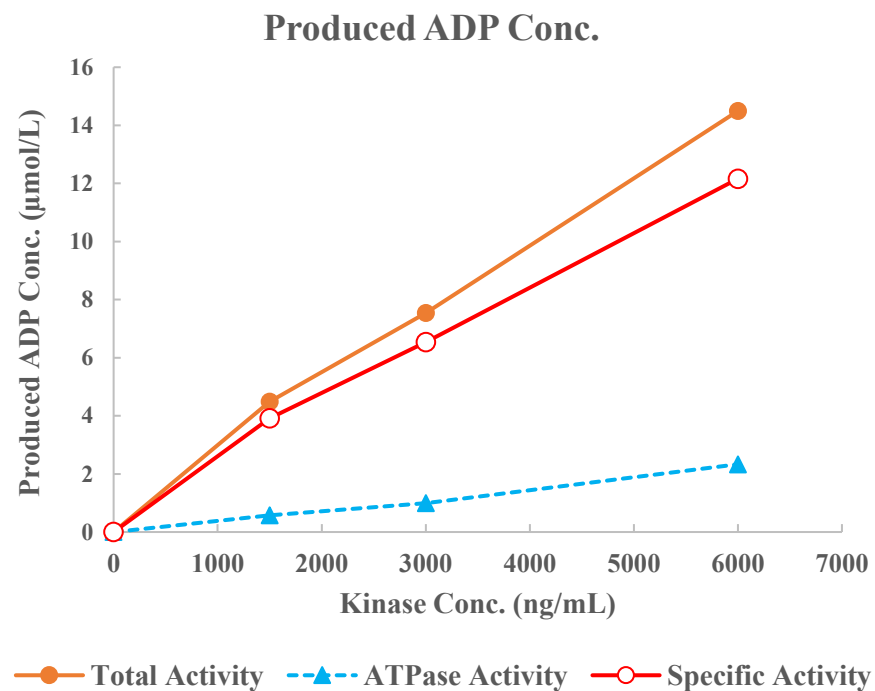
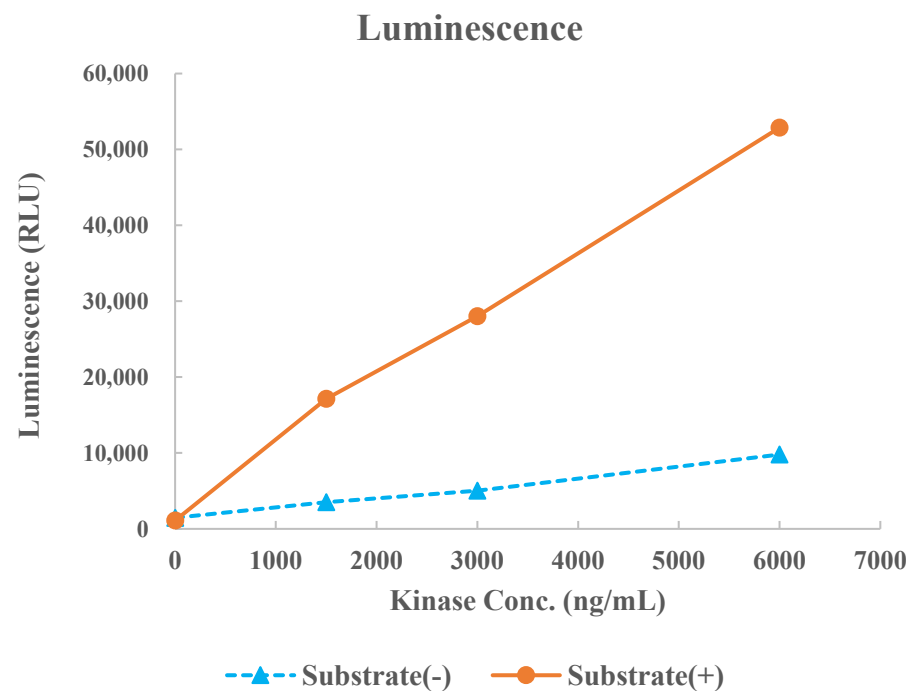


MSK2(RPS6KA4)



Kinase	MSK2(RPS6KA4), 01-148
Substrate	Biotin-Crosstide, BTSU19
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour