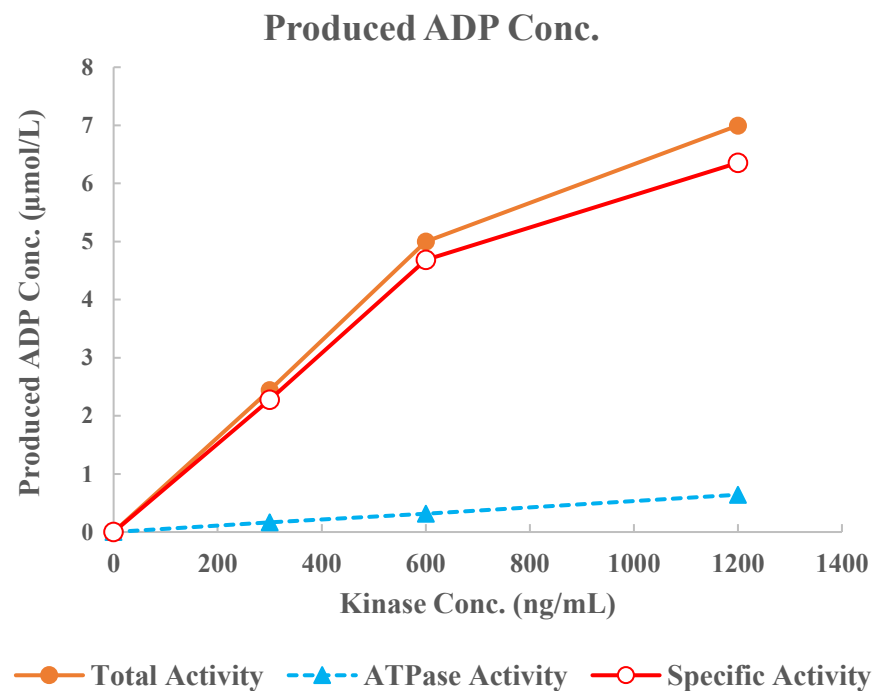
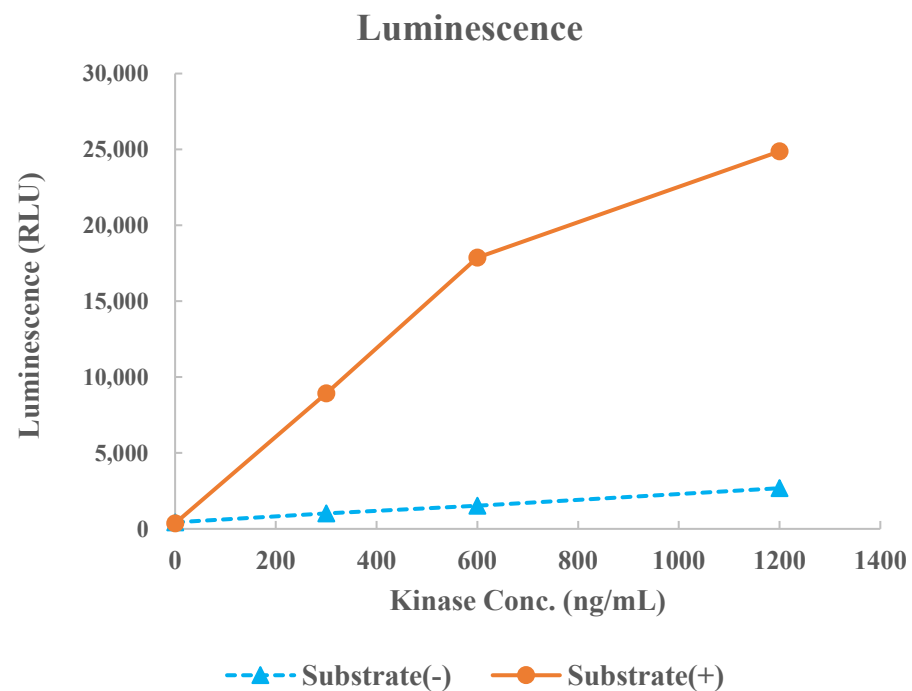


MSK1(RPS6KA5)



Kinase	MSK1(RPS6KA5), 01-147
Substrate	Biotin-Crosstide, BTSU19
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour