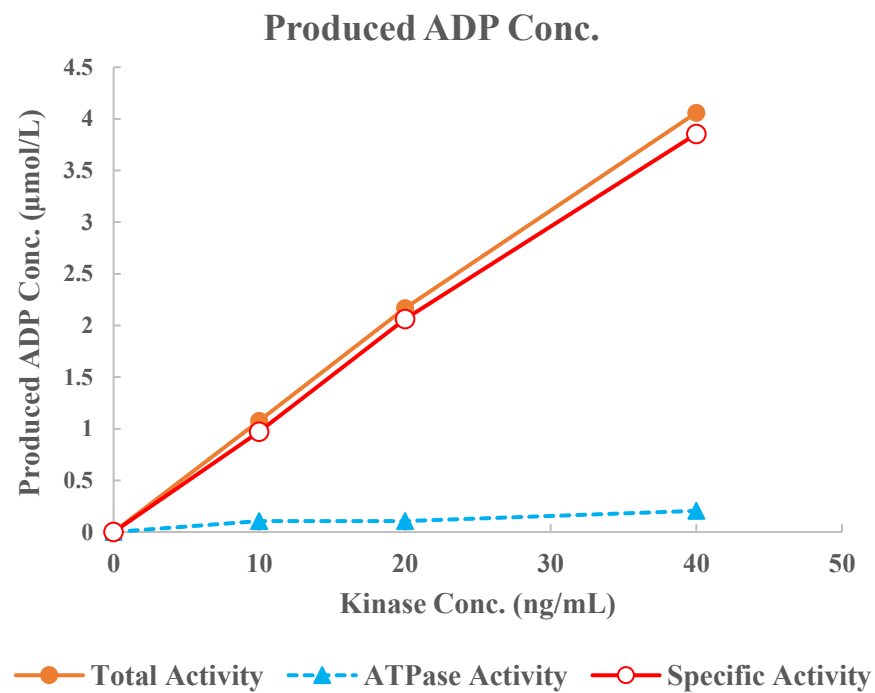
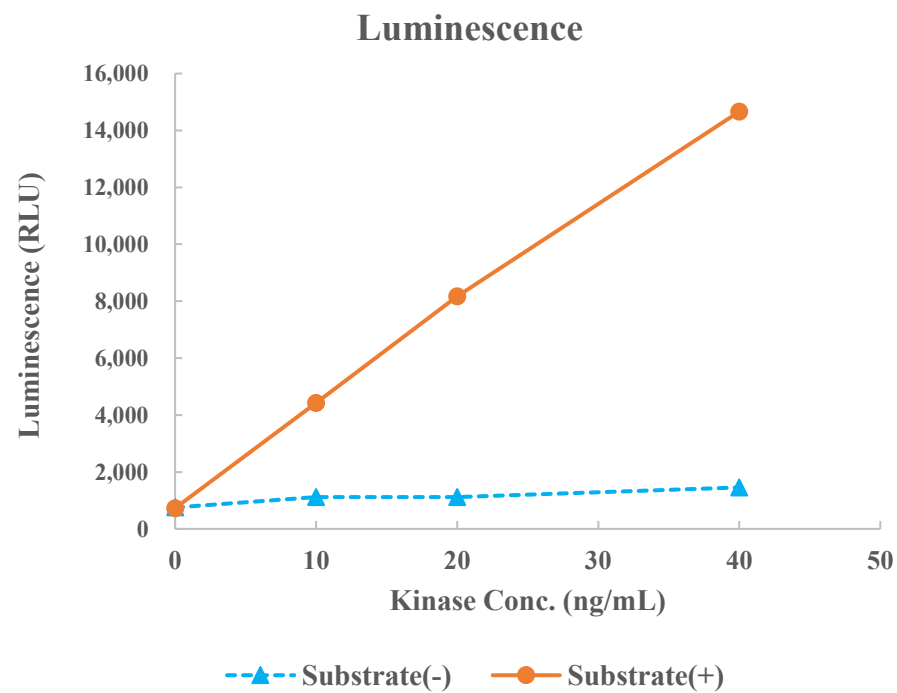


RSK4(RPS6KA6)



Kinase	RSK4(RPS6KA6), 01-152
Substrate	Biotin-S6K peptide, BTSU20
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour