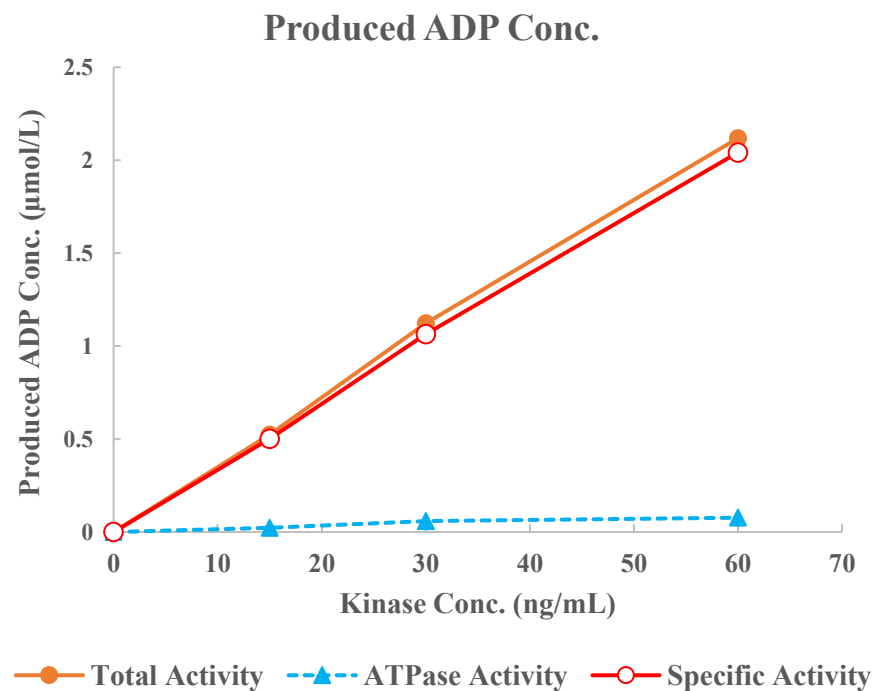
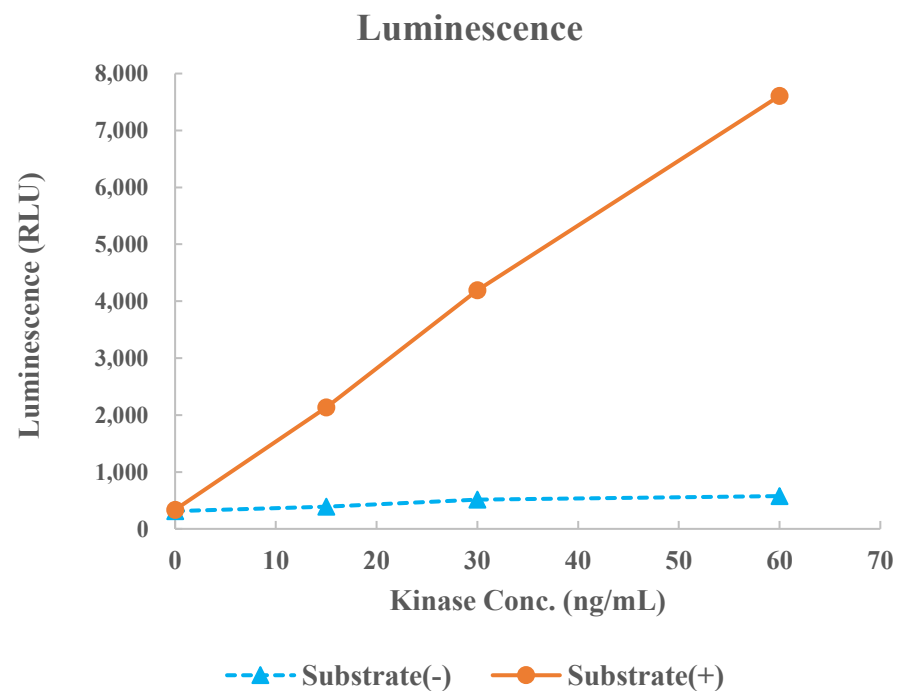


RSK2(RPS6KA3)



Kinase	RSK2(RPS6KA3), 01-150
Substrate	Biotin-S6K peptide, BTSU20
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour