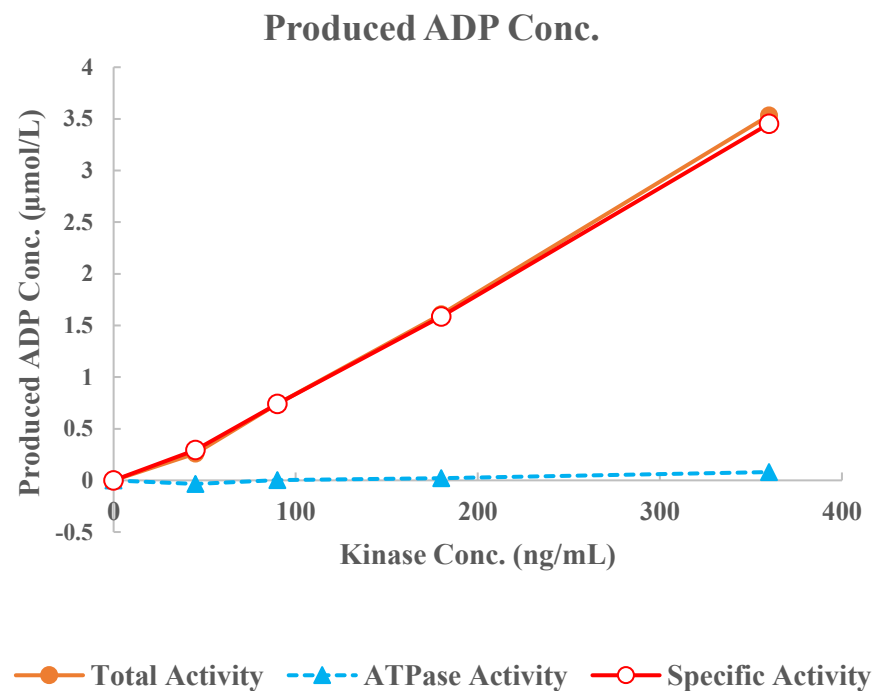
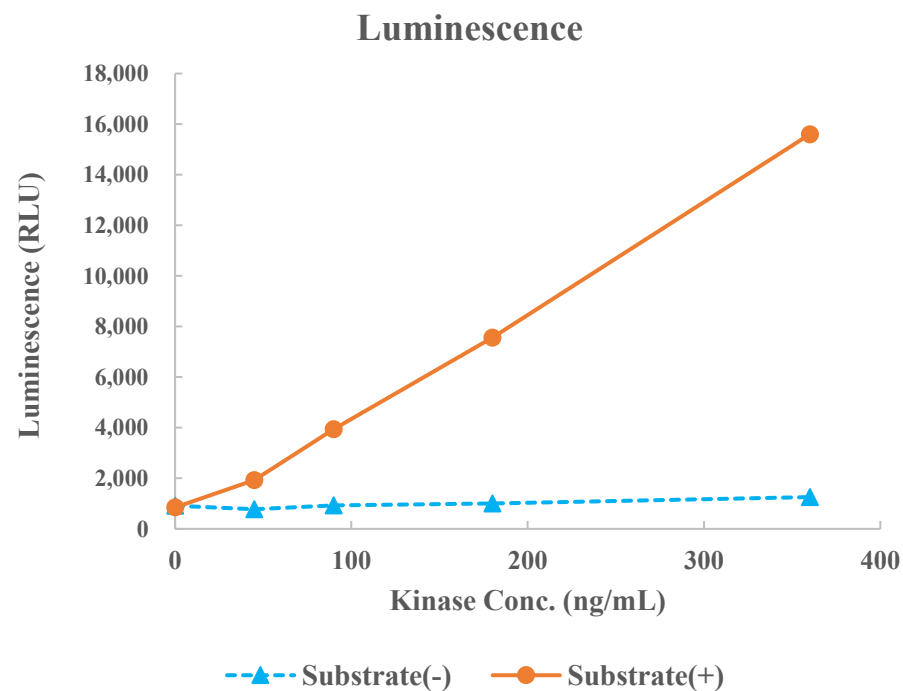


RSK1(RPS6KA1)



Kinase	RSK1(RPS6KA1), 01-149
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