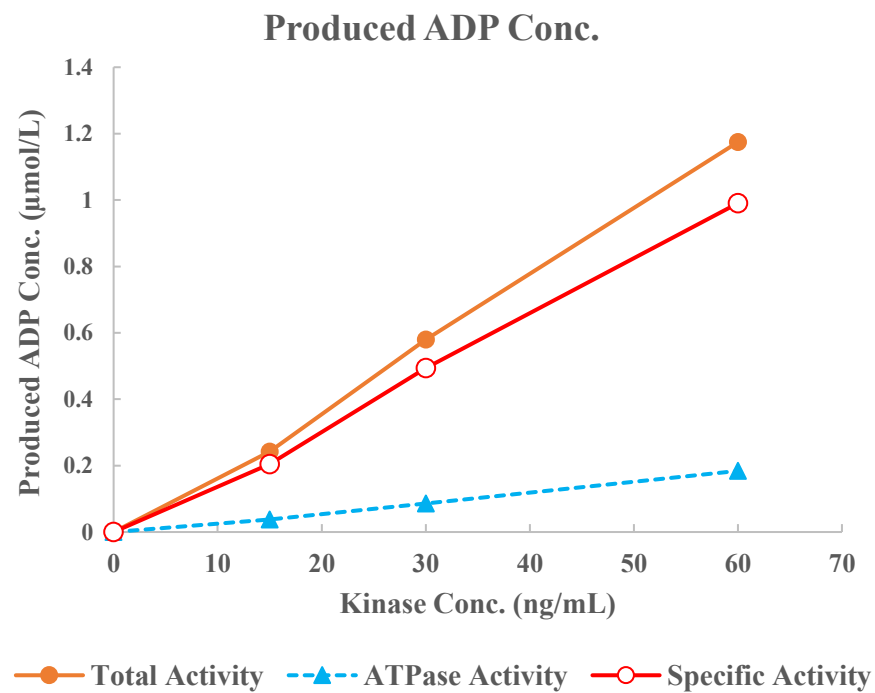
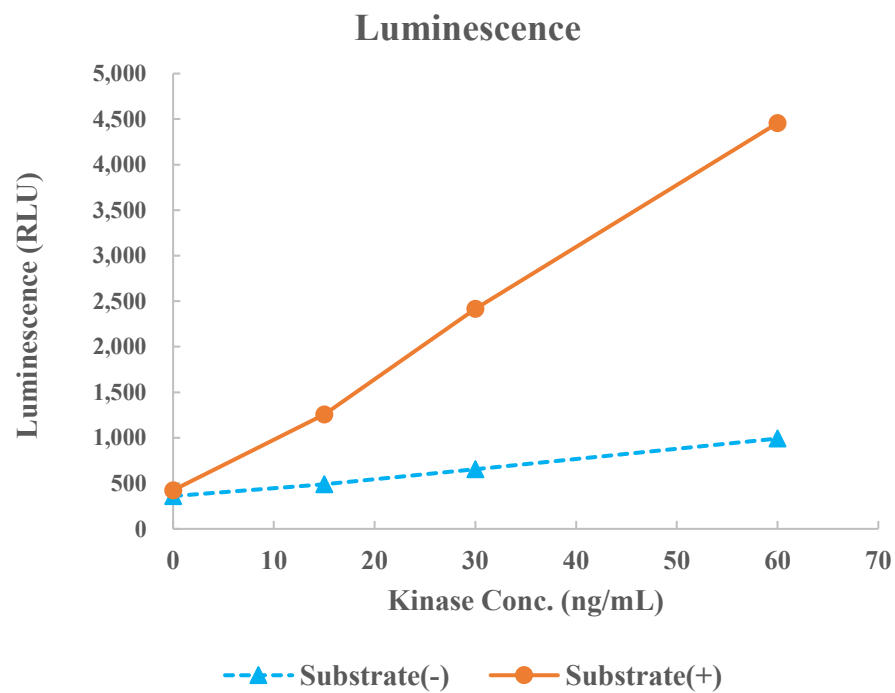


RSK3(RPS6KA2)



Kinase	RSK3(RPS6KA2), 01-151
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