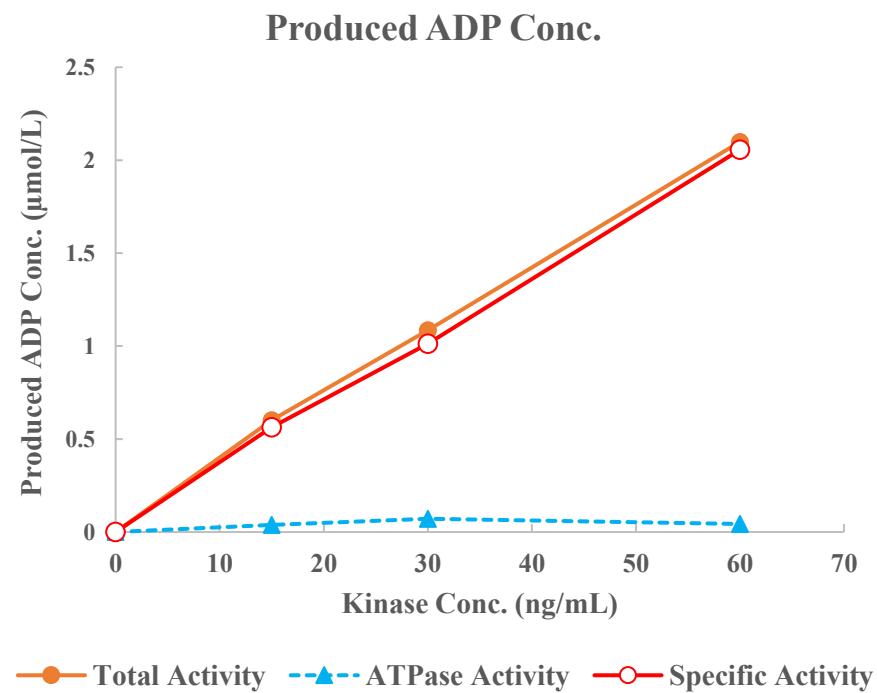
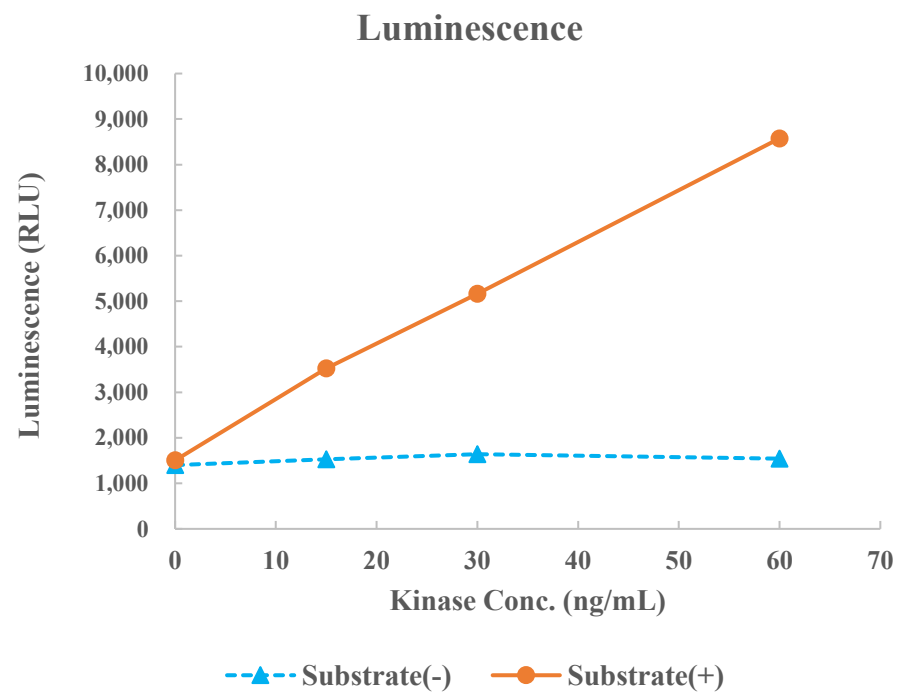


SGK2



Kinase	SGK2, 01-159
Substrate	Biotin-SGKtide, BTSU15
Substrate Conc.	10 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour