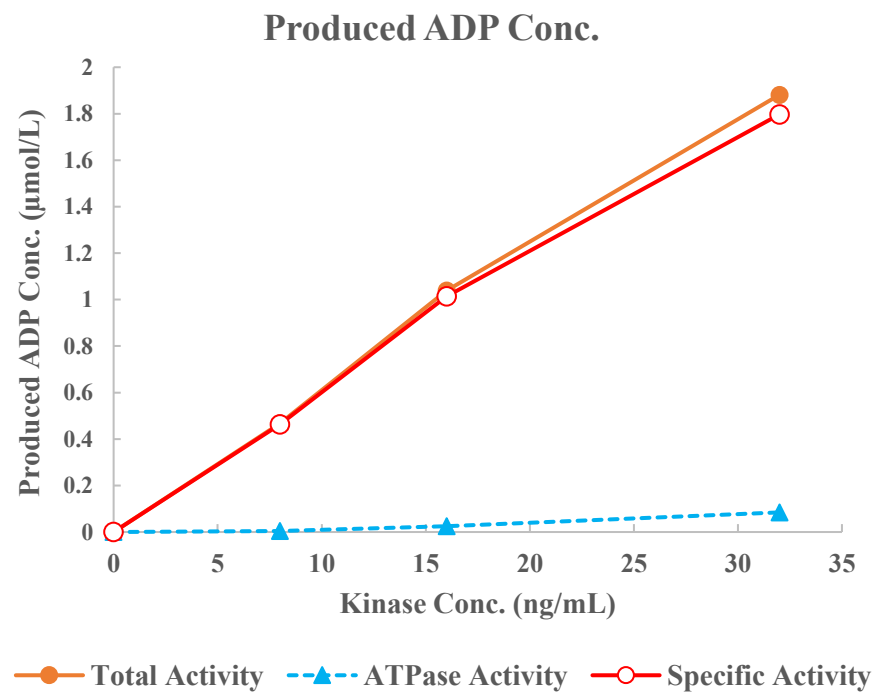
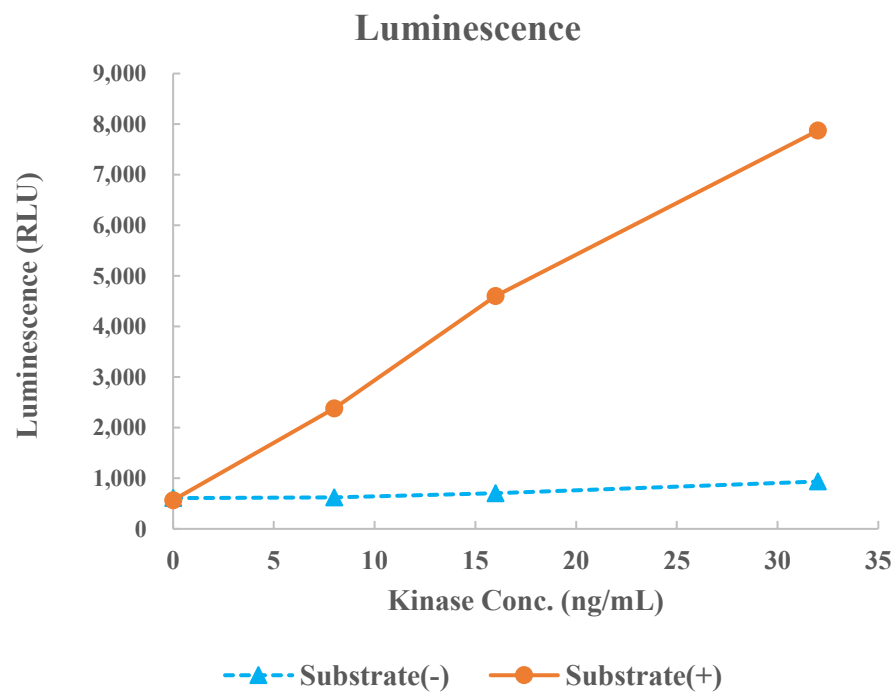


SGK3(SGKL)



Kinase	SGK3(SGKL), 01-160
Substrate	Biotin-SGKtide, BTSU15
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour