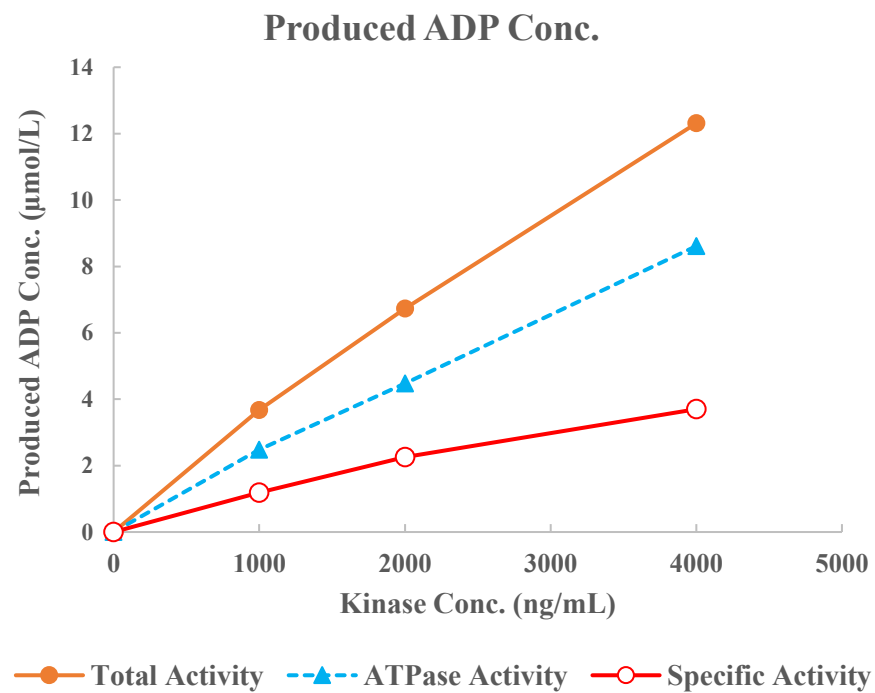
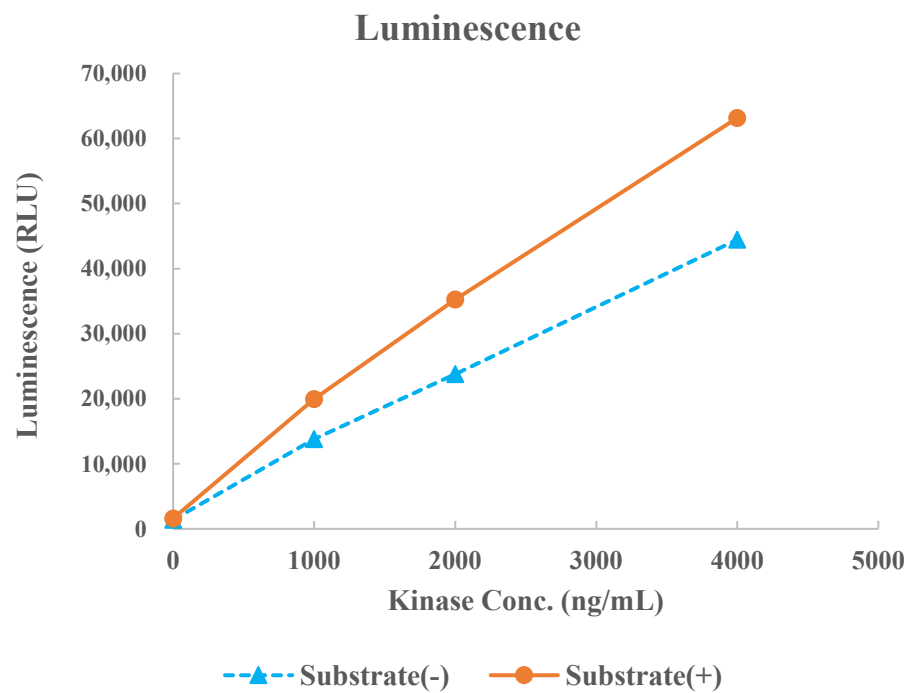


WNK2



Kinase	WNK2, 05-180
Substrate	Biotin-SPAKtide, BTSU27
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg, Mn
Metal Conc.	5, 3 mmol/L
Additives	-
Reaction Time	1 hour