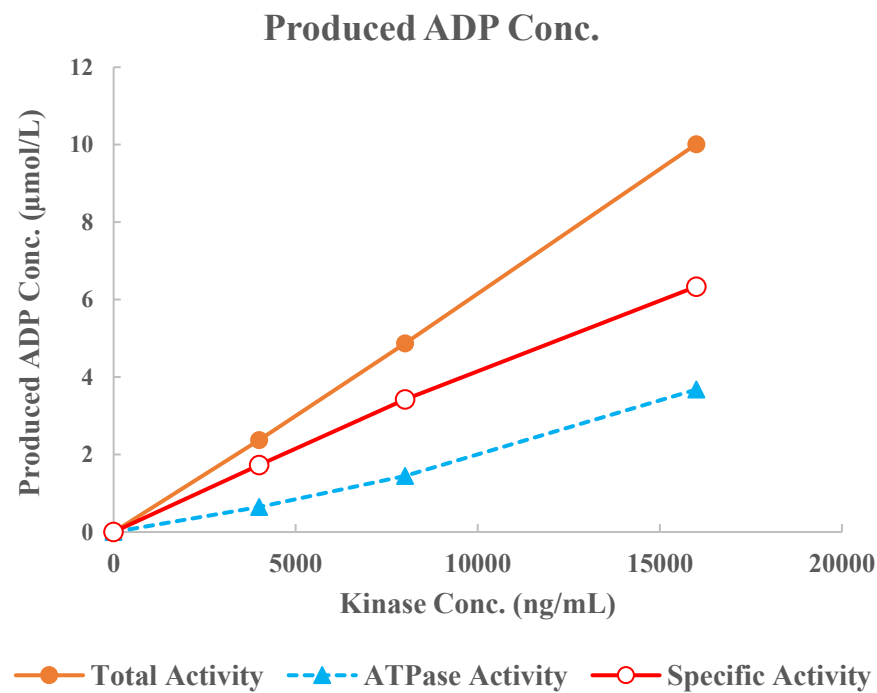
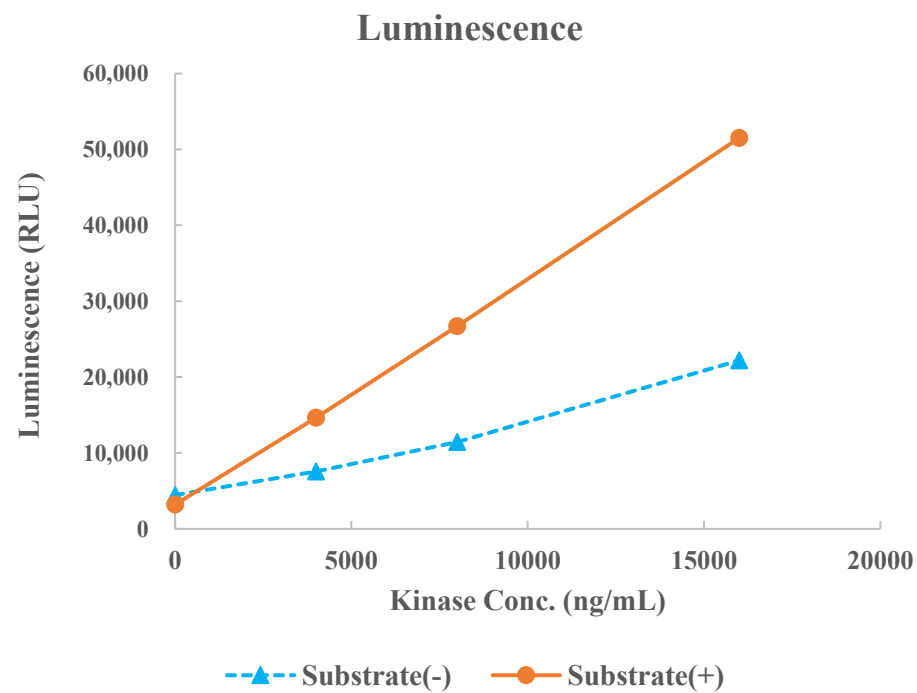


BTN-WNK4



Kinase	BTN-WNK4, 05-482-20N
Substrate	Biotin-SPAKtide, BTSU27
Substrate Conc.	50 μmol/L
ATP Conc.	150 μmol/L
Metal	Mg, Mn
Metal Conc.	5, 3 mmol/L
Additives	-
Reaction Time	1 hour