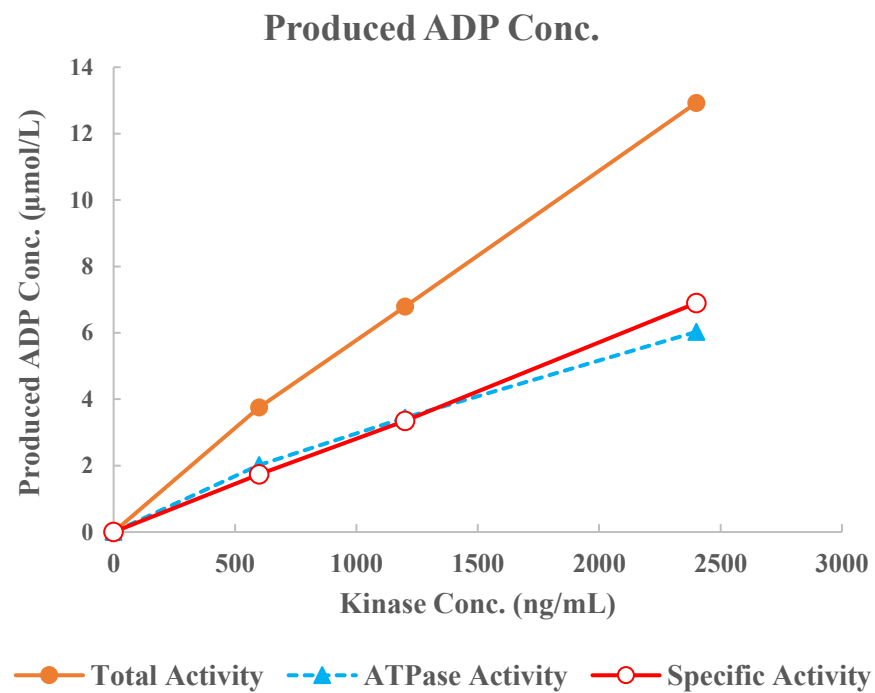
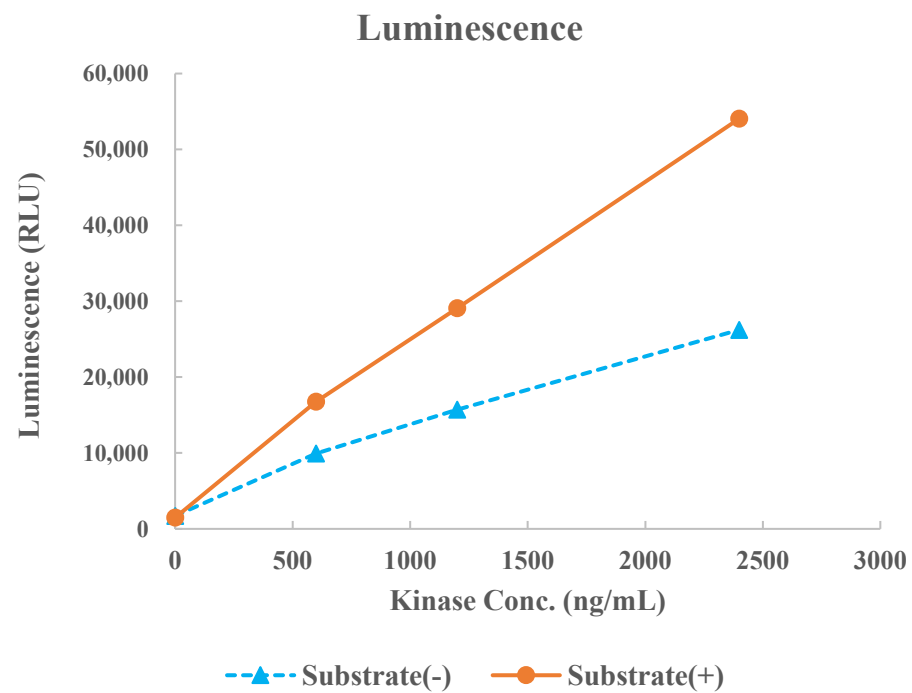


TNK1



Kinase	TNK1, 08-104
Substrate	Biotin-CSKtide, BTSU02
Substrate Conc.	50 μmol/L
ATP Conc.	75 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour