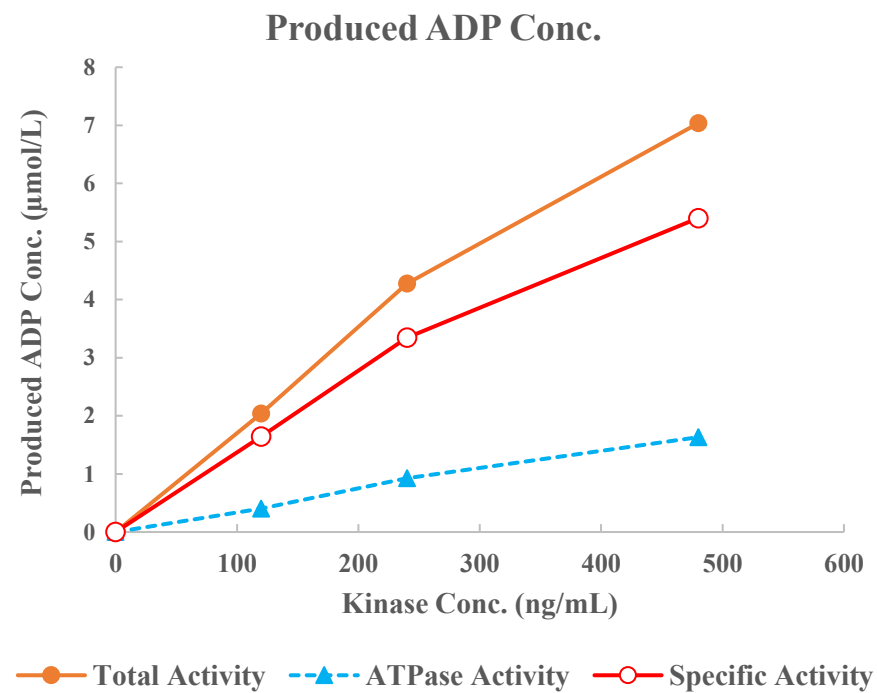
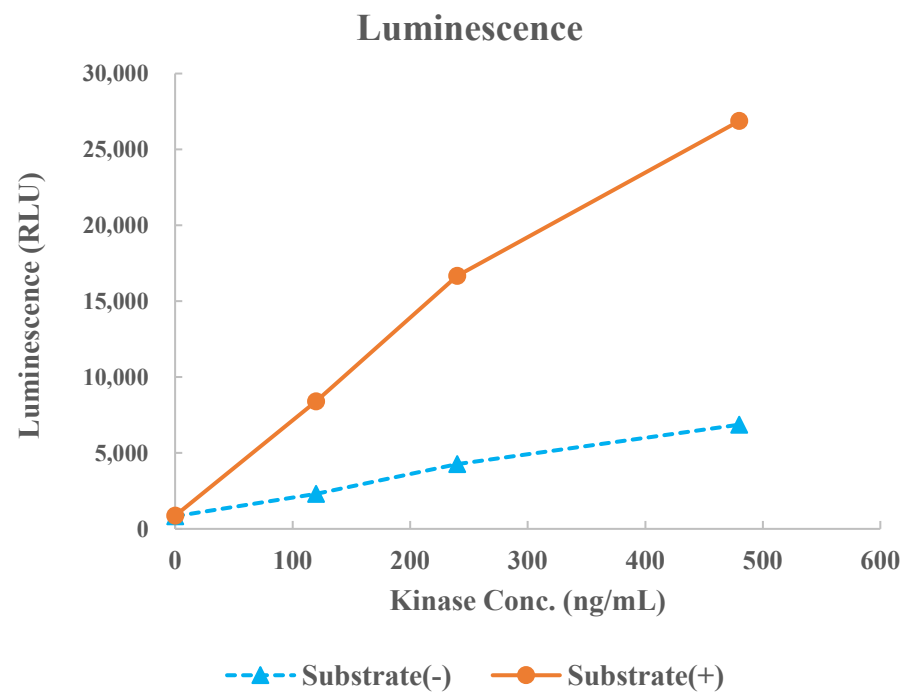


TBK1



Kinase	TBK1, 05-115
Substrate	Biotin-CKtide, BTSU12
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour