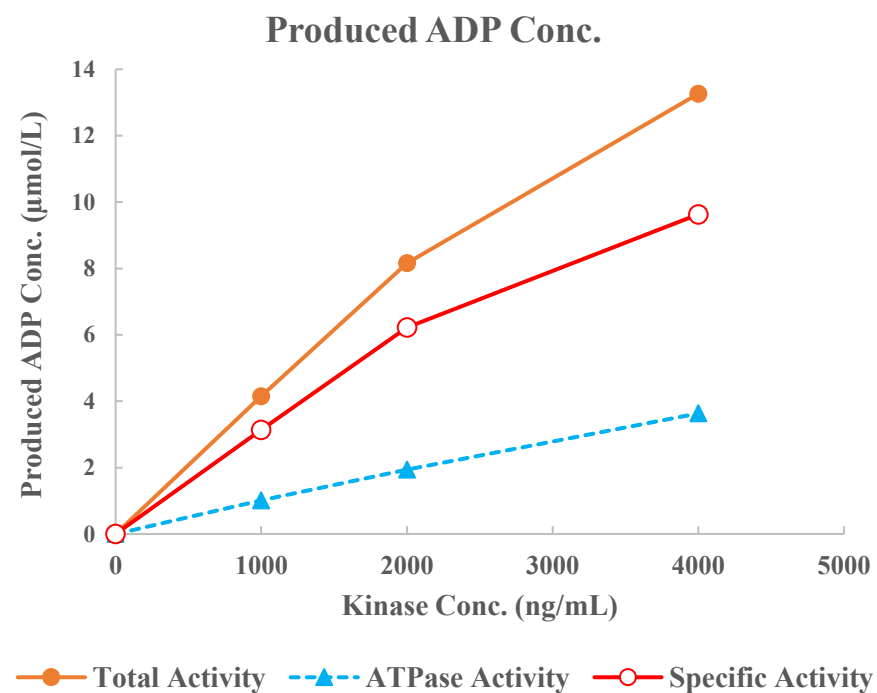
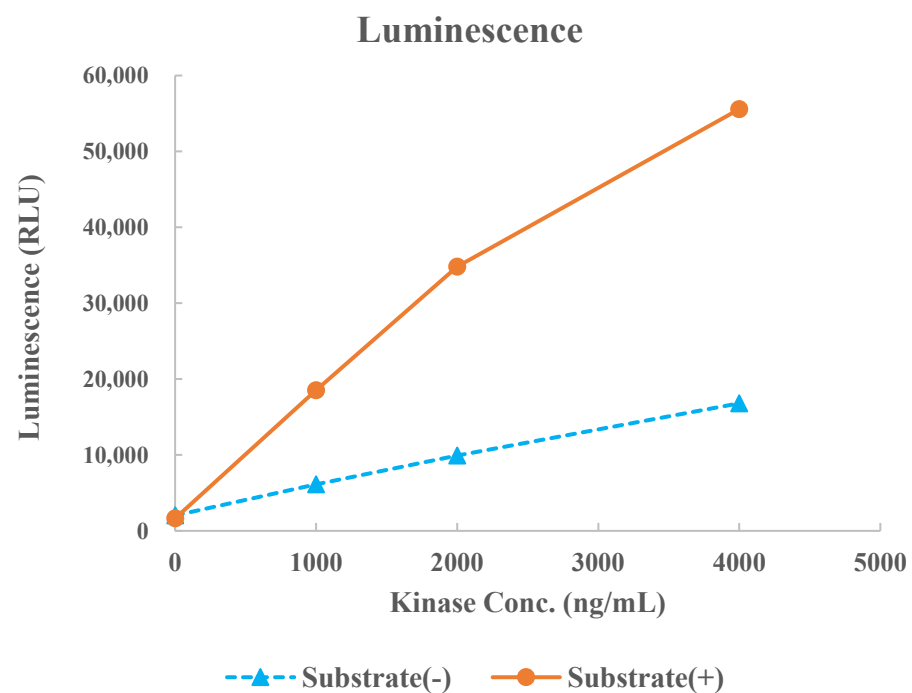


NEK6



Kinase	NEK6, 05-130
Substrate	Biotin-CDK7 peptide, BTSU11
Substrate Conc.	50 μmol/L
ATP Conc.	75 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour