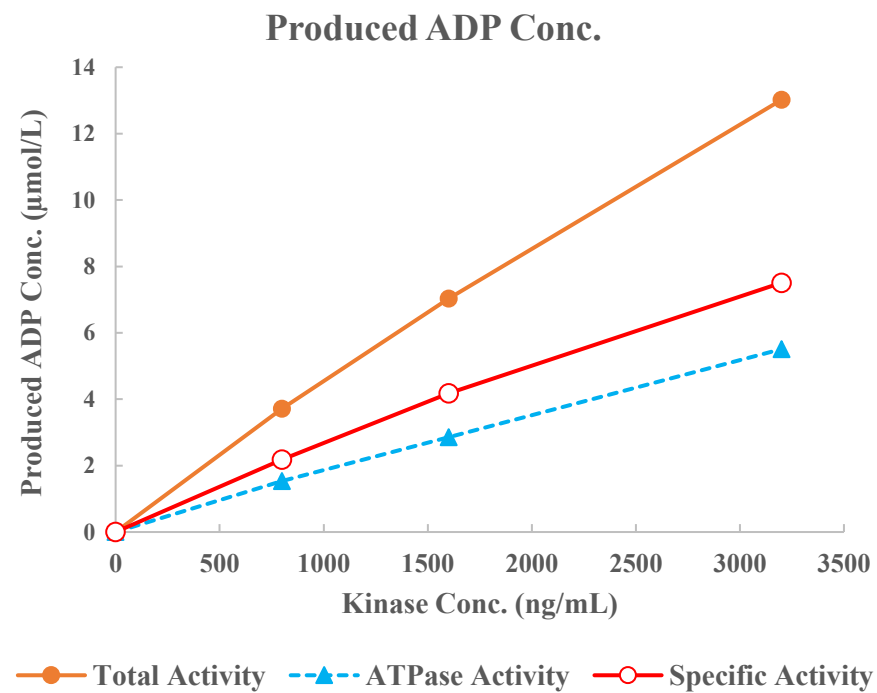
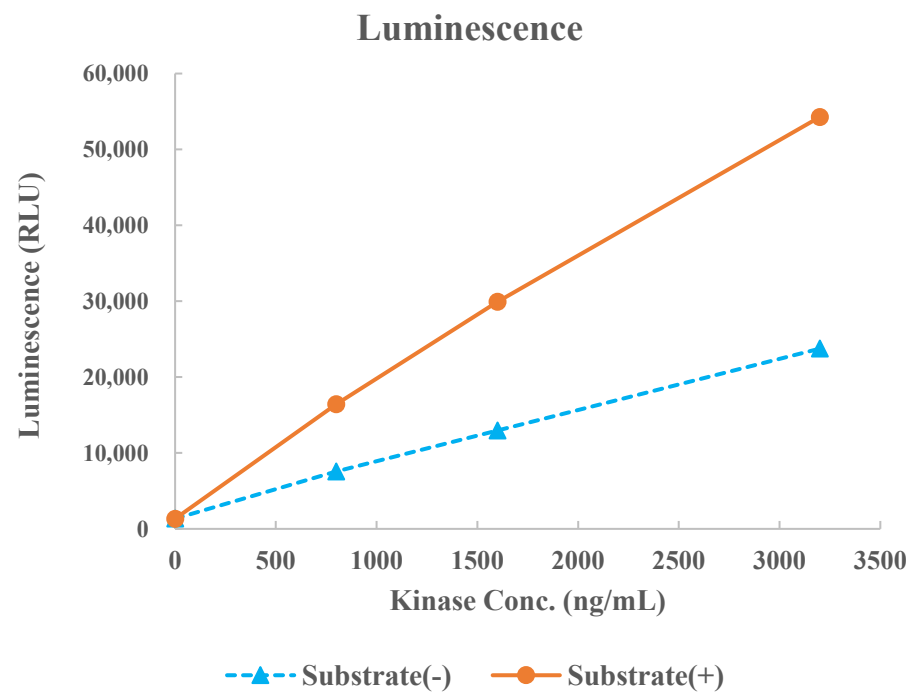


NEK7



Kinase	NEK7, 05-131
Substrate	Biotin-CDK7 peptide, BTSU11
Substrate Conc.	50 μmol/L
ATP Conc.	50 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour