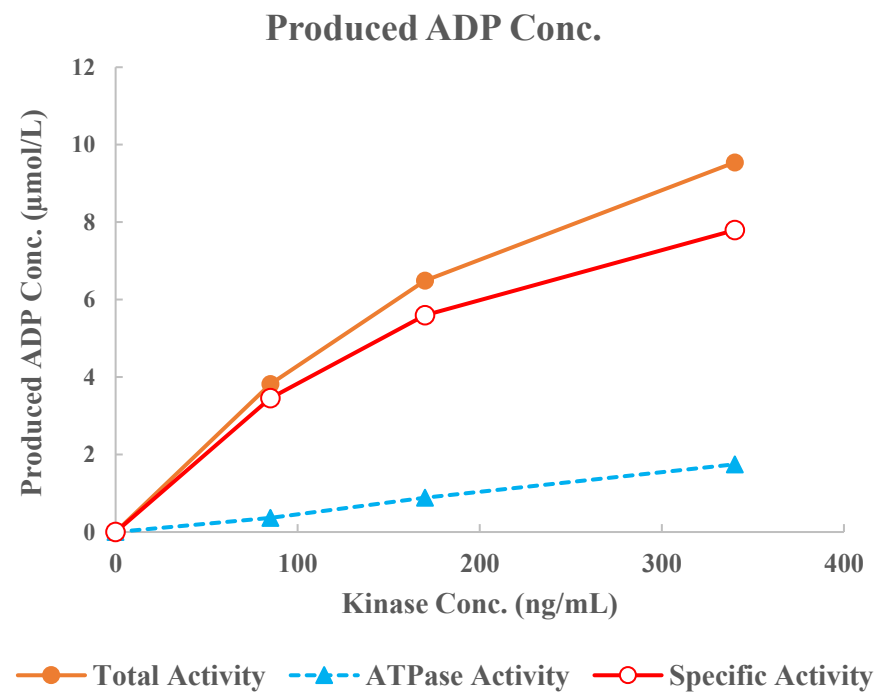
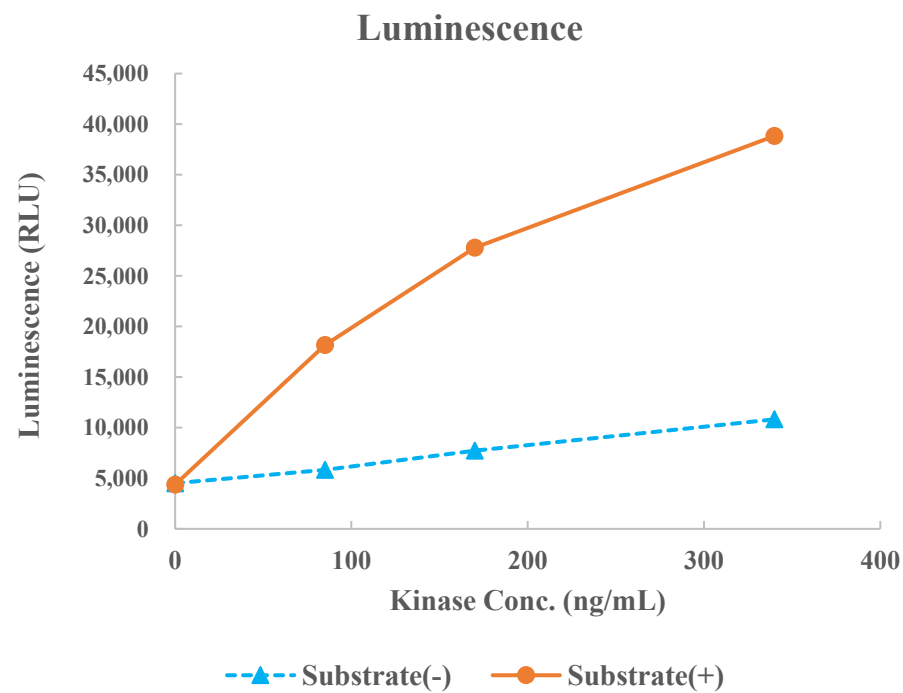


NEK9



Kinase	NEK9, 05-133
Substrate	Biotin-CDK7 peptide, BTSU11
Substrate Conc.	10 μmol/L
ATP Conc.	200 μmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 hour