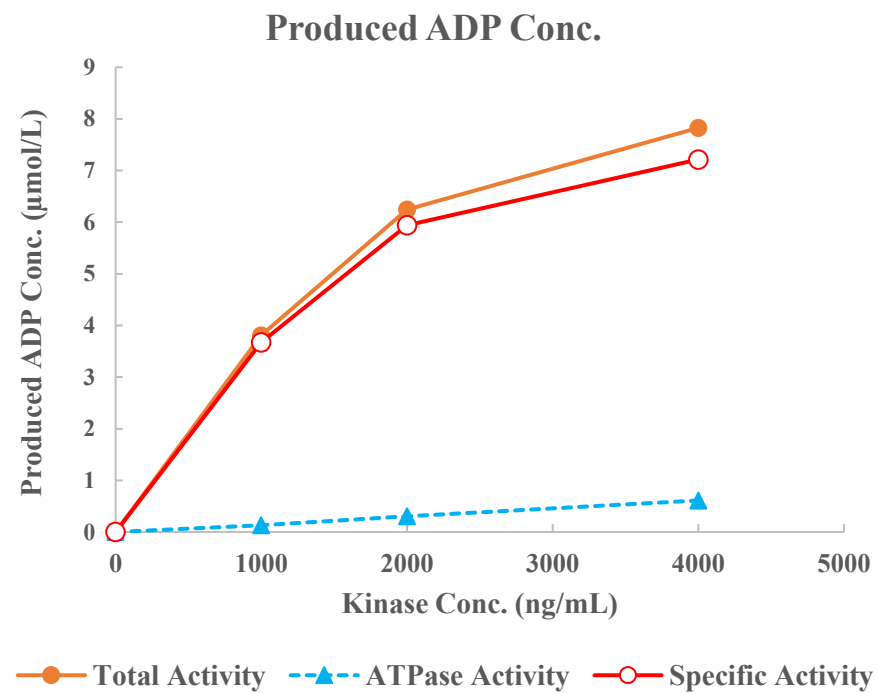
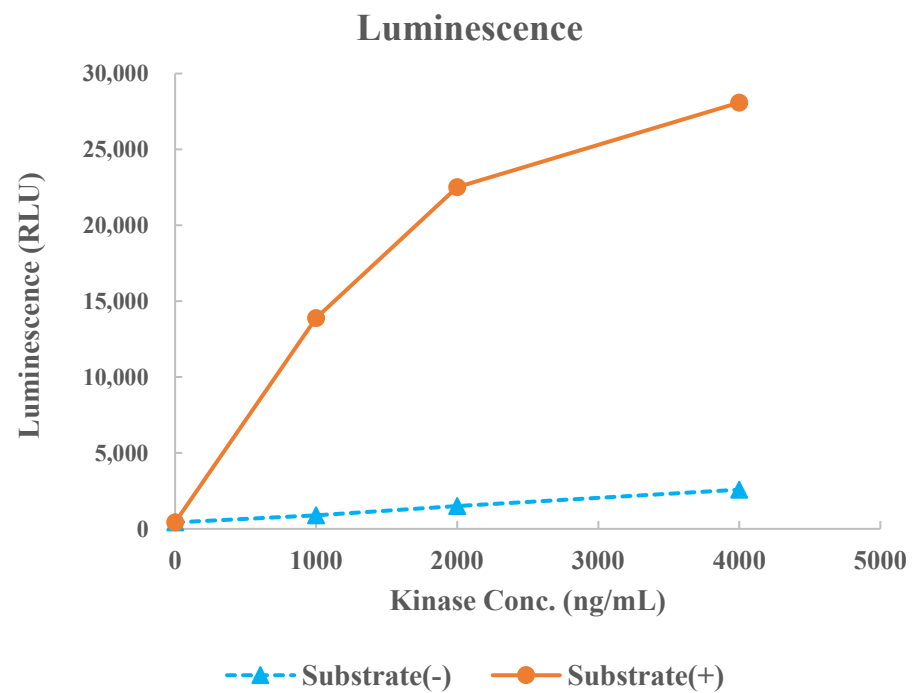


EEF2K



Kinase	EEF2K, 10-113
Substrate	Biotin-EEF2Ktide, BTSU40
Substrate Conc.	10 μmol/L
ATP Conc.	10 μmol/L
Metal	Mg, Ca
Metal Conc.	5, 1 mmol/L
Additives	10 μg/mL Calmodulin
Reaction Time	1 hour