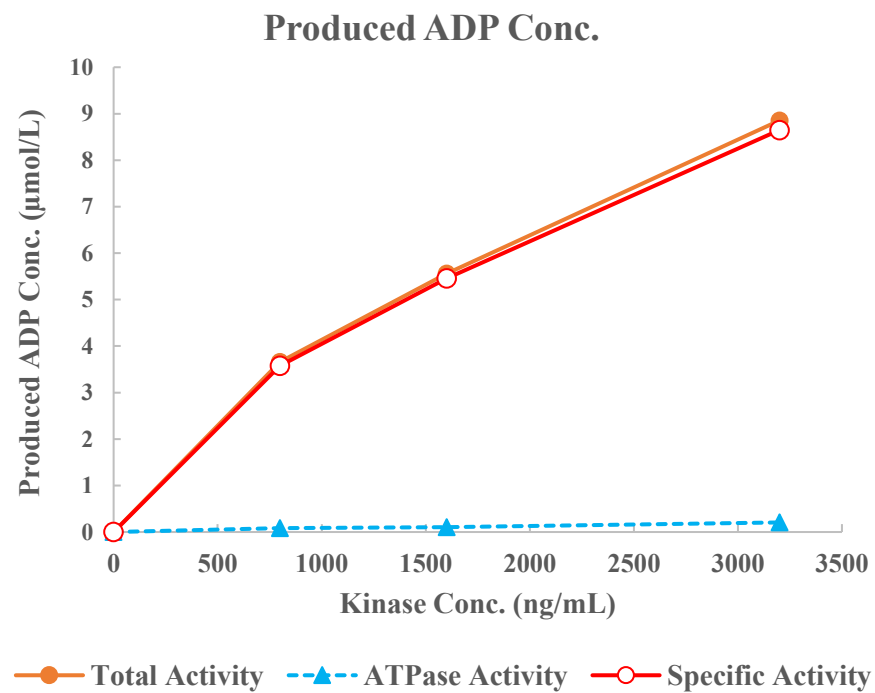
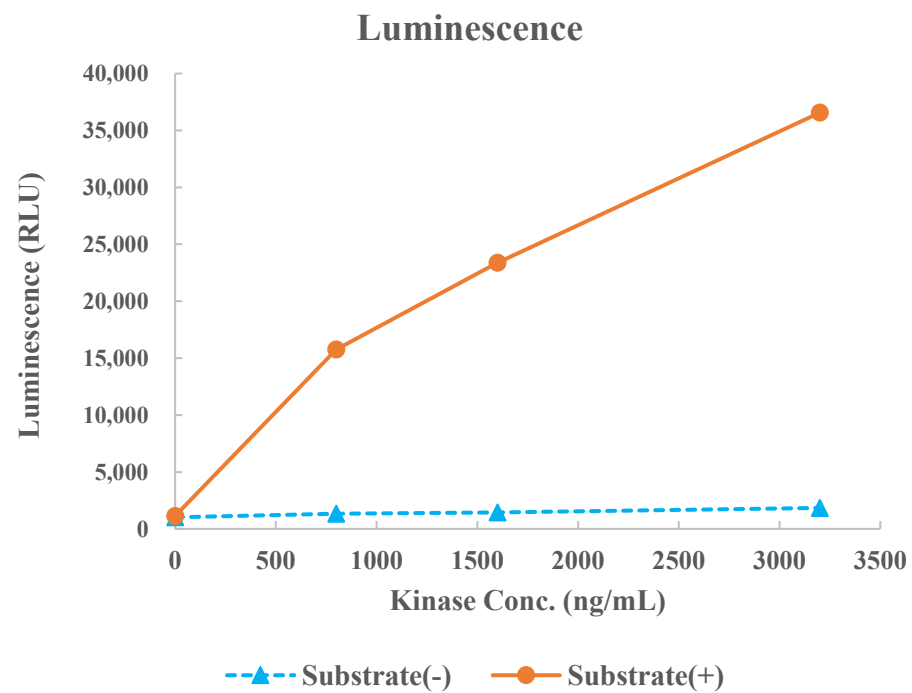


CaMK4



Kinase	CaMK4, 02-108
Substrate	Biotin-GS peptide, BTSU05
Substrate Conc.	10 μmol/L
ATP Conc.	25 μmol/L
Metal	Mg, Ca
Metal Conc.	5, 1 mmol/L
Additives	10 μg/mL Calmodulin
Reaction Time	1 hour