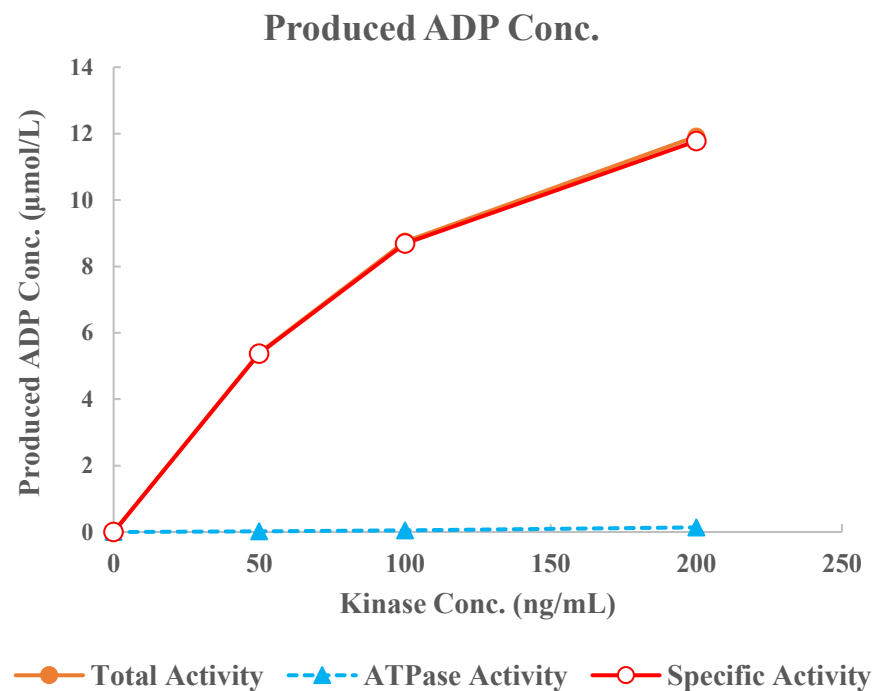
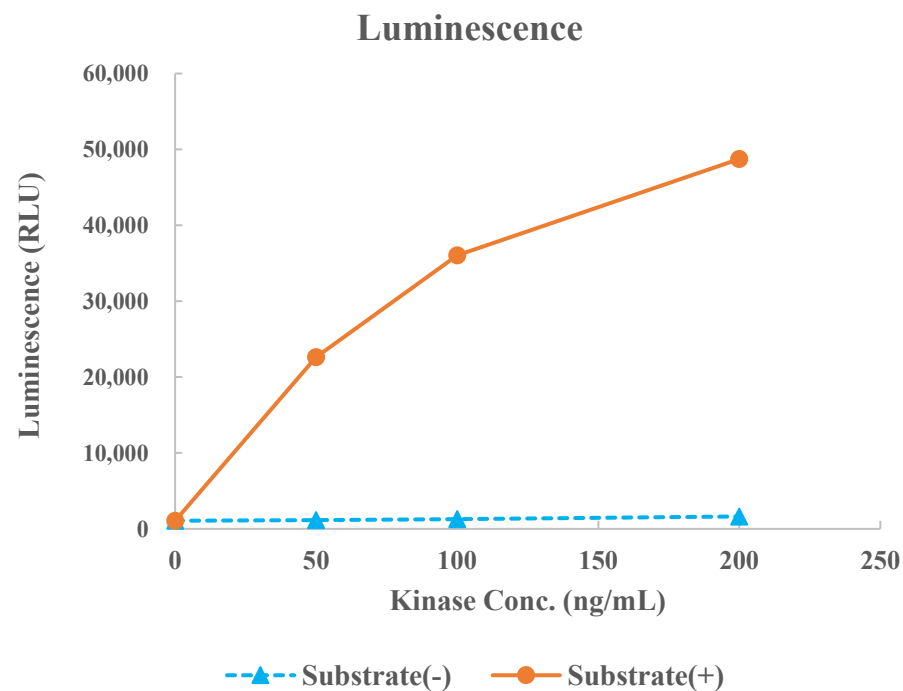


CaMK2 β (CAMK2B)



Kinase
 Substrate
 Substrate Conc.
 ATP Conc.
 Metal
 Metal Conc.
 Additives
 Reaction Time

CaMK2 β (CAMK2B), 02-110
 Biotin-GS peptide, BTSU05
 10 μ mol/L
 25 μ mol/L
 Mg, Ca
 5, 1 mmol/L
 10 μ g/mL Calmodulin
 1 hour