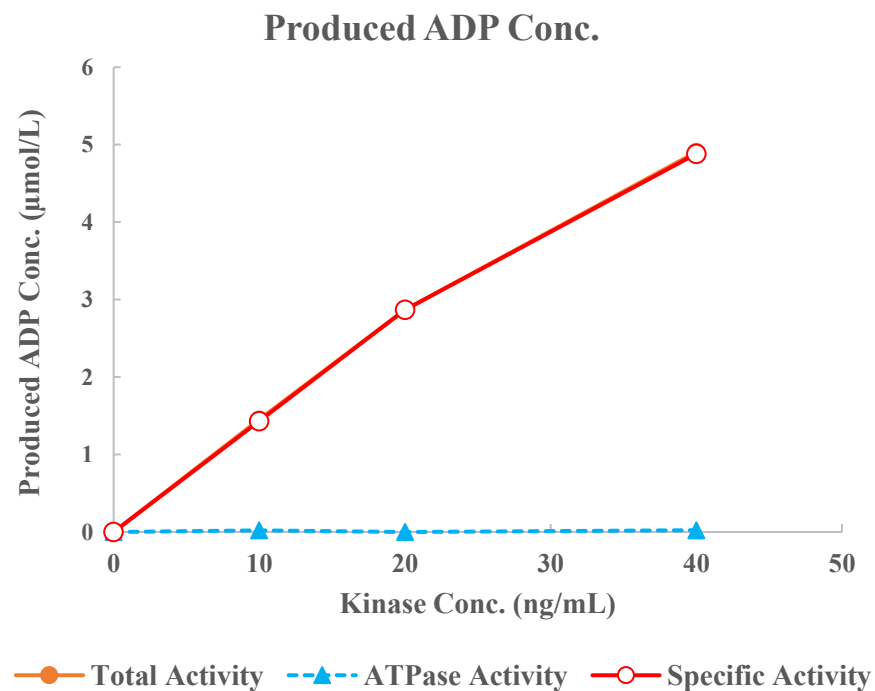
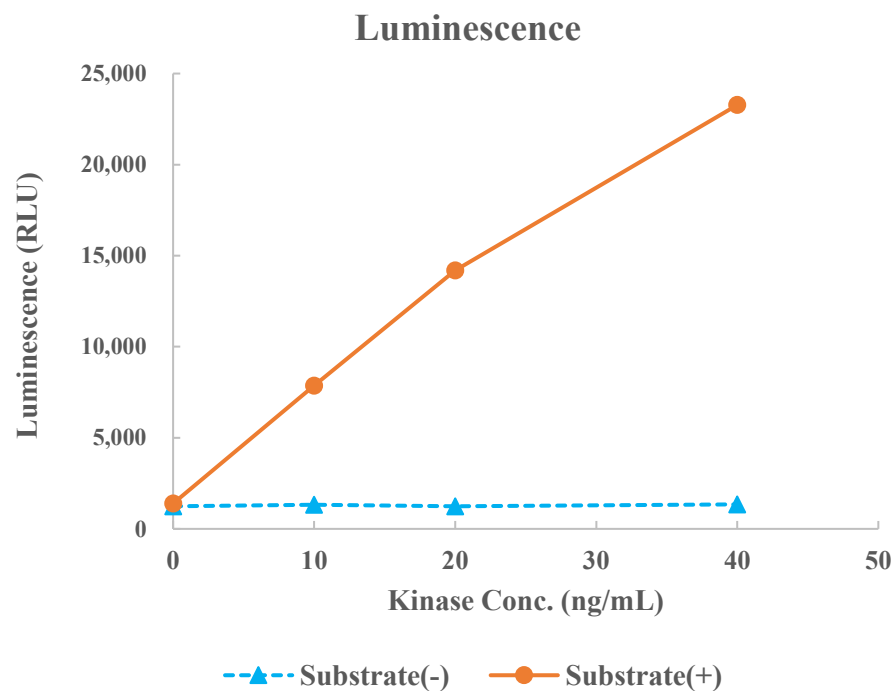


PKCθ(PRKCQ)



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

PKCθ(PRKCQ), 01-140
 Biotin-PKC peptide, BTSU10
 10 μmol/L
 25 μmol/L
 Mg
 5 mmol/L
 5 μg/mL Diacyl Glycerol, 50 μg/mL Phosphatidylserine
 1 hour