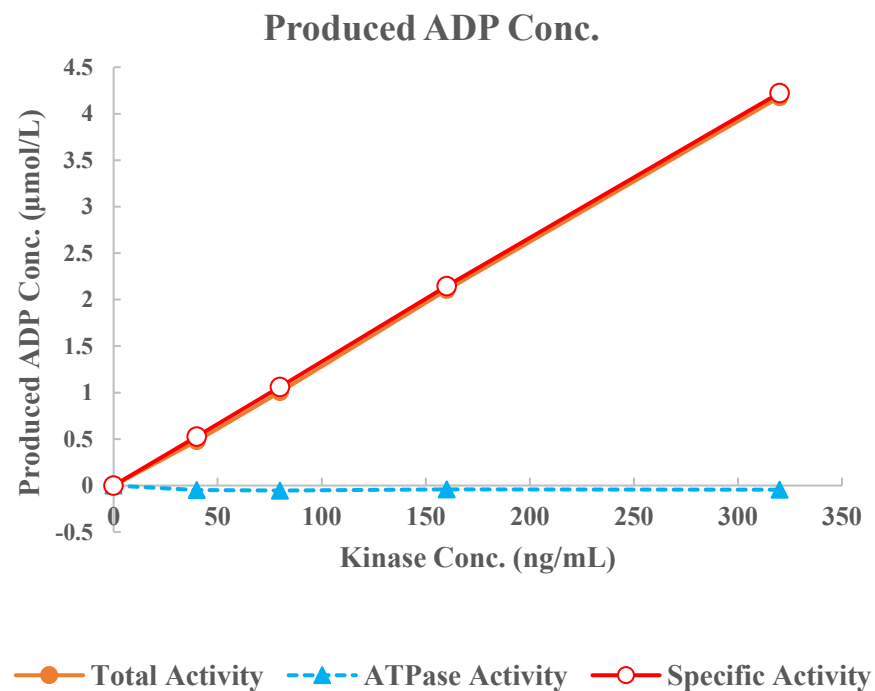
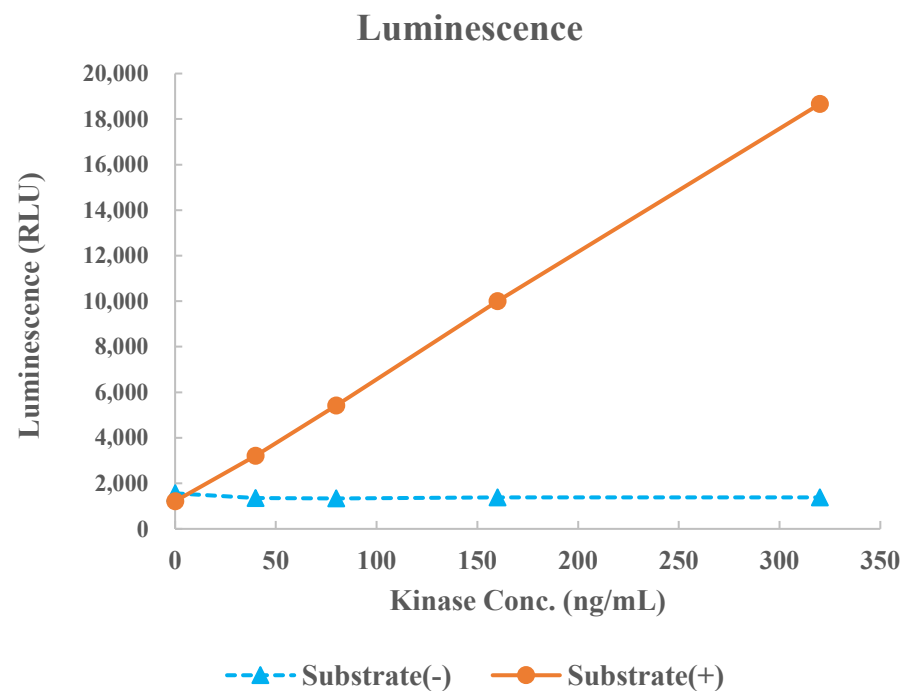


PKC η (PRKCH)



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

PKC η (PRKCH), 01-138
 Biotin-PKC peptide, BTSU10
 10 μ mol/L
 50 μ mol/L
 Mg
 5 mmol/L
 5 μ g/mL Diacyl Glycerol, 50 μ g/mL Phosphatidylserine
 1 hour