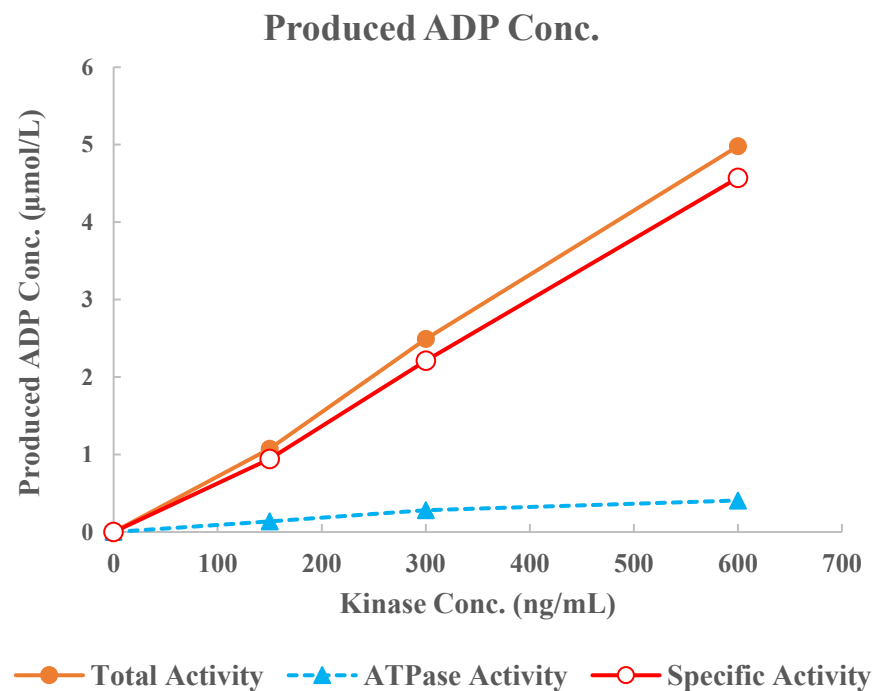
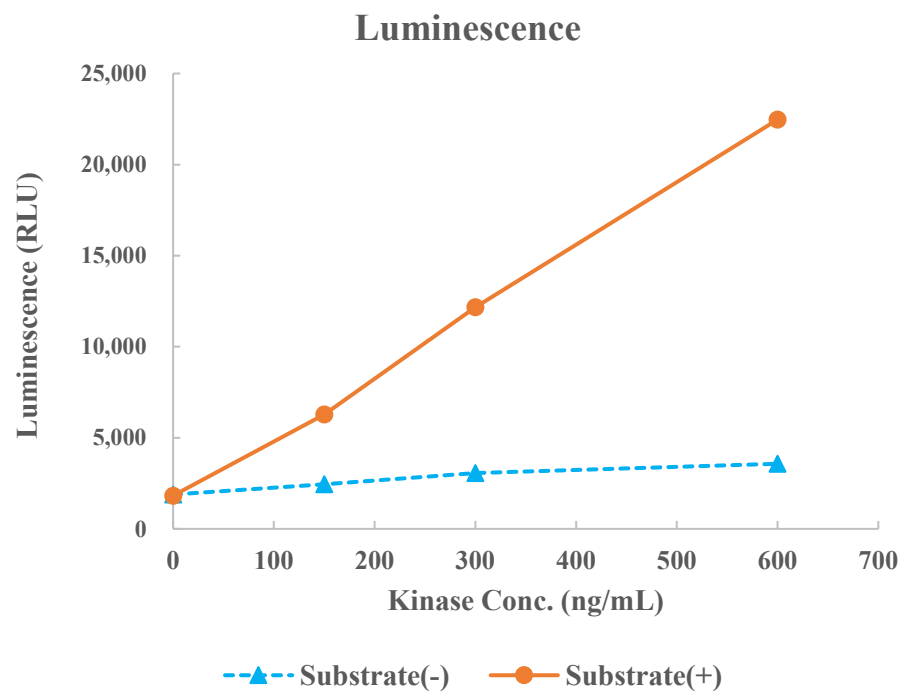


PDGFR α (PDGFRA) [V561D]



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

PDGFR α (PDGFRA) [V561D], 08-507
 Biotin-CSKtide, BTSU02
 10 μ mol/L
 50 μ mol/L
 Mg
 5 mmol/L
 -
 1 hour