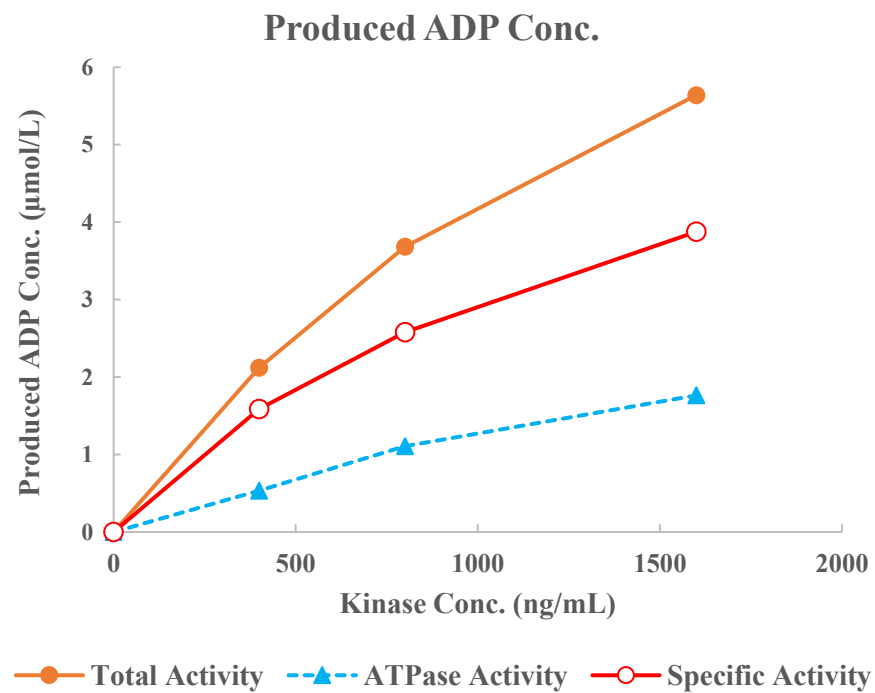
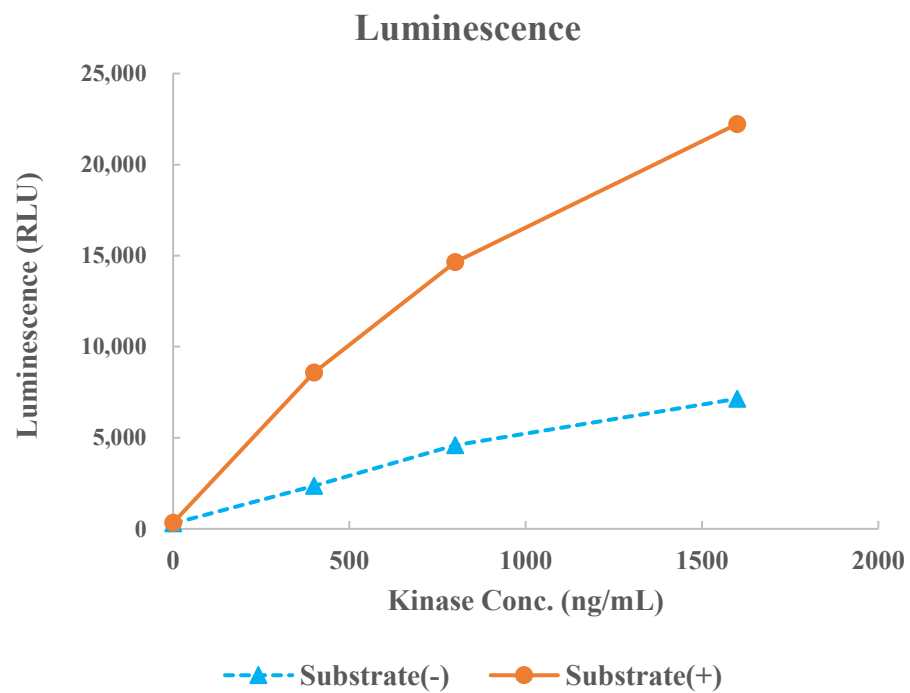


## EGFR [T790M/C797S/L858R]



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

EGFR [T790M/C797S/L858R], 08-559  
 Biotin-Srcptide, BTSU01  
 10 μmol/L  
 10 μmol/L  
 Mg, Mn  
 5, 1 mmol/L  
 -  
 1 hour