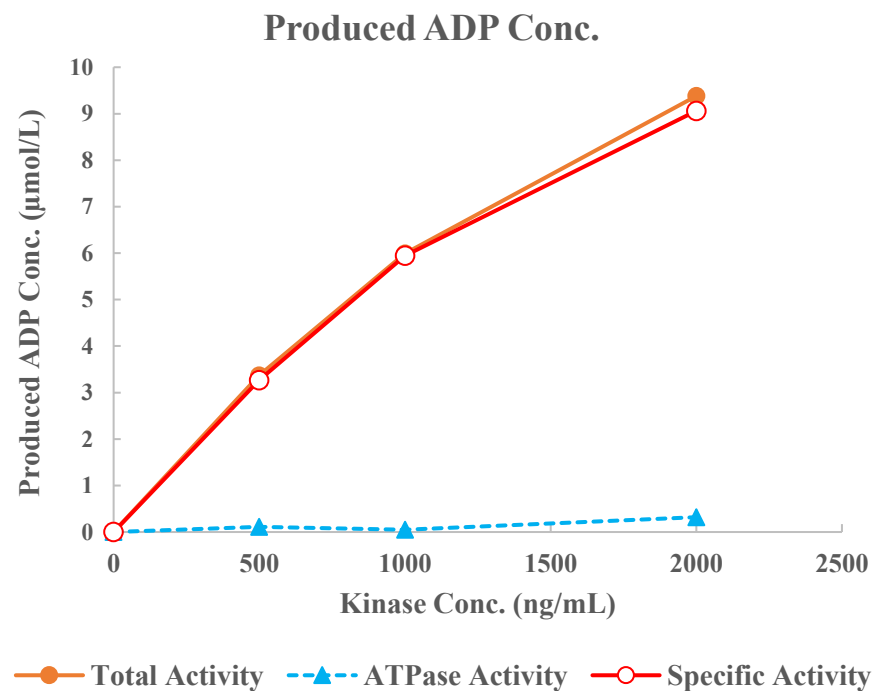
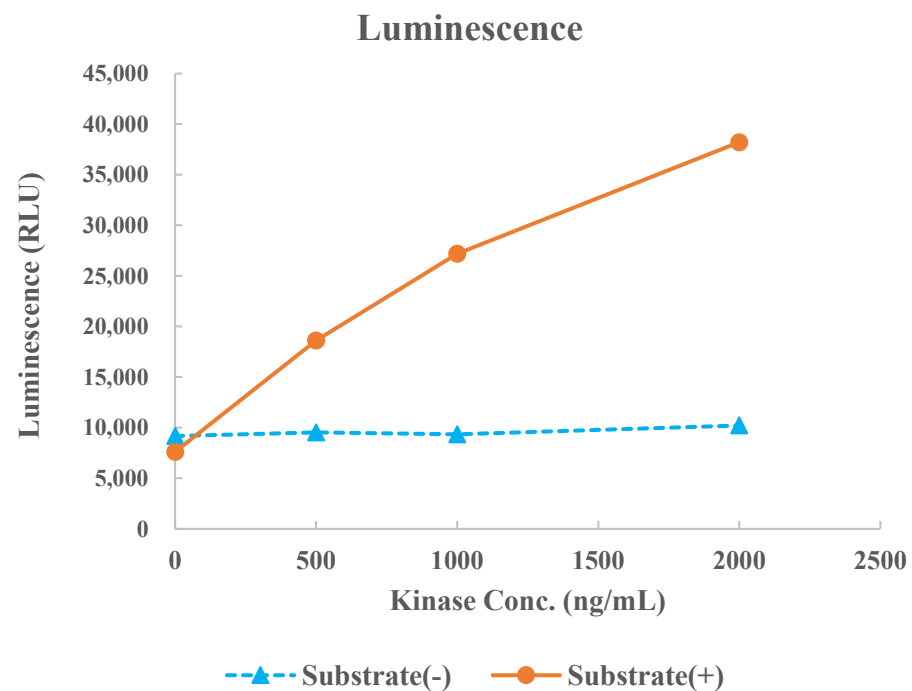


## KIT [V560G]



|                 |                                |
|-----------------|--------------------------------|
| Kinase          | KIT [V560G], 08-504            |
| Substrate       | Biotin-Srcptide, BTSU01        |
| Substrate Conc. | 10 μmol/L                      |
| ATP Conc.       | 250 μmol/L                     |
| Metal           | Mg                             |
| Metal Conc.     | 5 mmol/L                       |
| Additives       | 25 μmol/L Sodium orthovanadate |
| Reaction Time   | 1 hour                         |