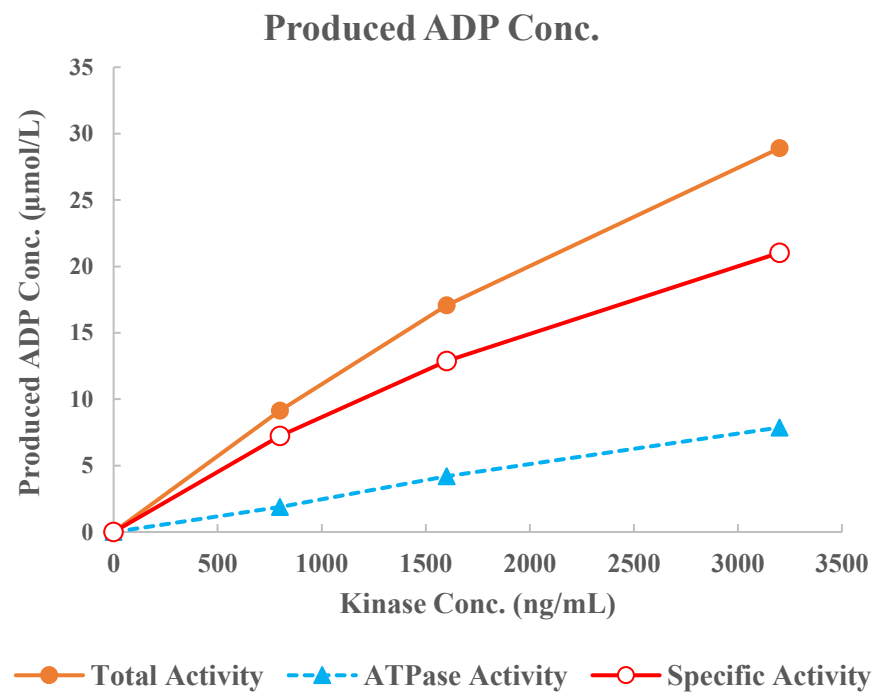
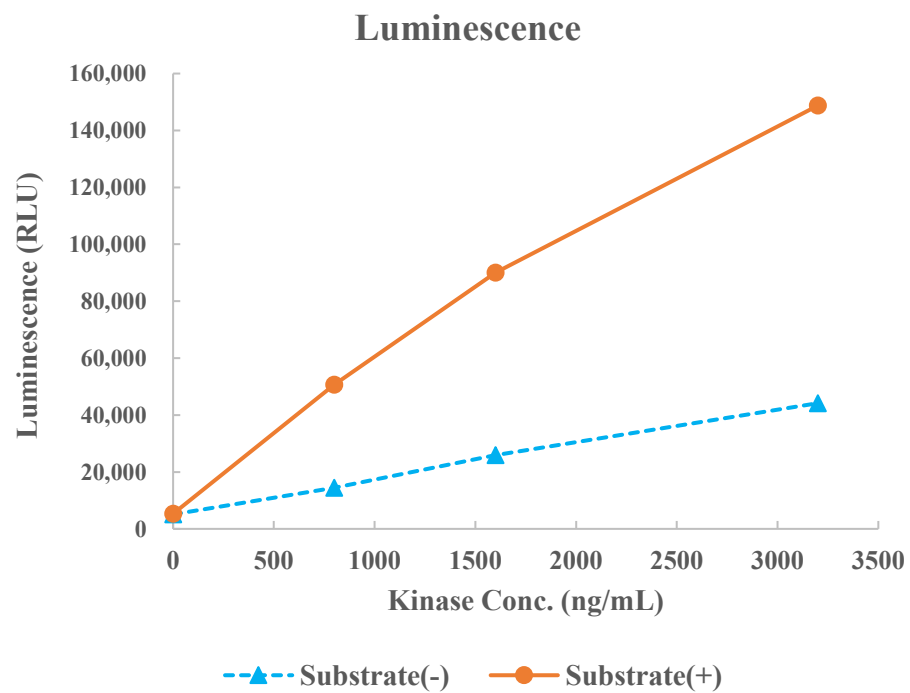


## ACK(TNK2)



|                 |                             |
|-----------------|-----------------------------|
| Kinase          | ACK(TNK2), 08-196           |
| Substrate       | Biotin-WASP peptide, BTSU39 |
| Substrate Conc. | 50 μmol/L                   |
| ATP Conc.       | 100 μmol/L                  |
| Metal           | Mg                          |
| Metal Conc.     | 5 mmol/L                    |
| Additives       | -                           |
| Reaction Time   | 1 hour                      |