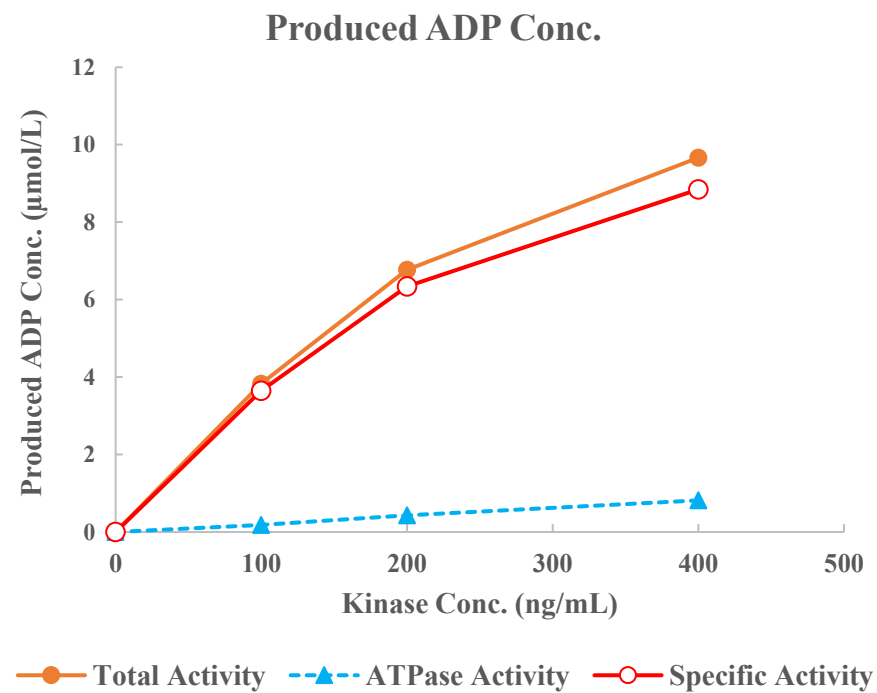
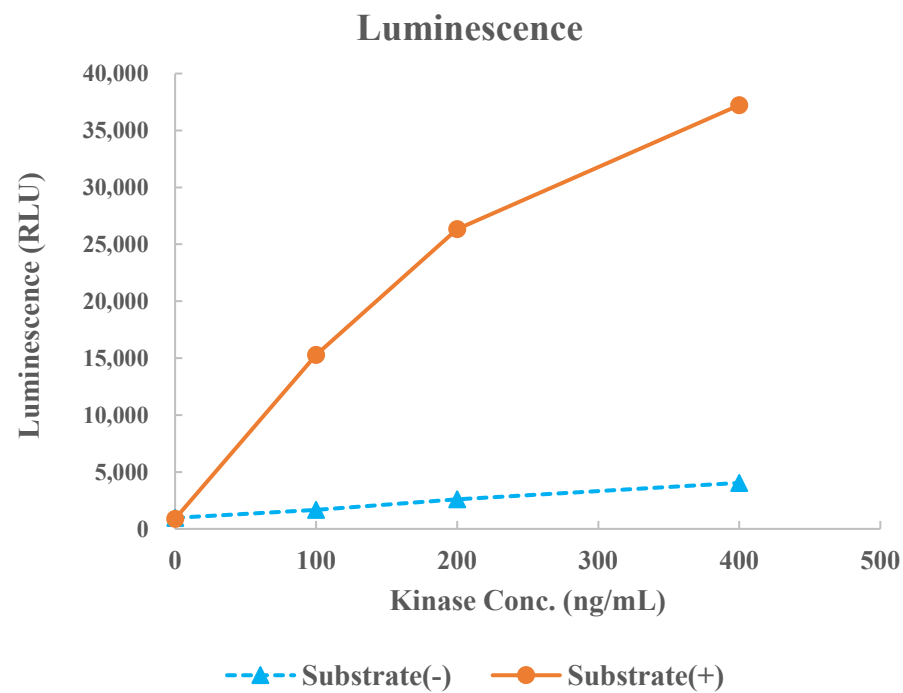


## DYRK1A



|                 |                         |
|-----------------|-------------------------|
| Kinase          | DYRK1A, 04-130          |
| Substrate       | Biotin-DYRKtide, BTSU06 |
| Substrate Conc. | 10 μmol/L               |
| ATP Conc.       | 25 μmol/L               |
| Metal           | Mg                      |
| Metal Conc.     | 5 mmol/L                |
| Additives       | -                       |
| Reaction Time   | 1 hour                  |