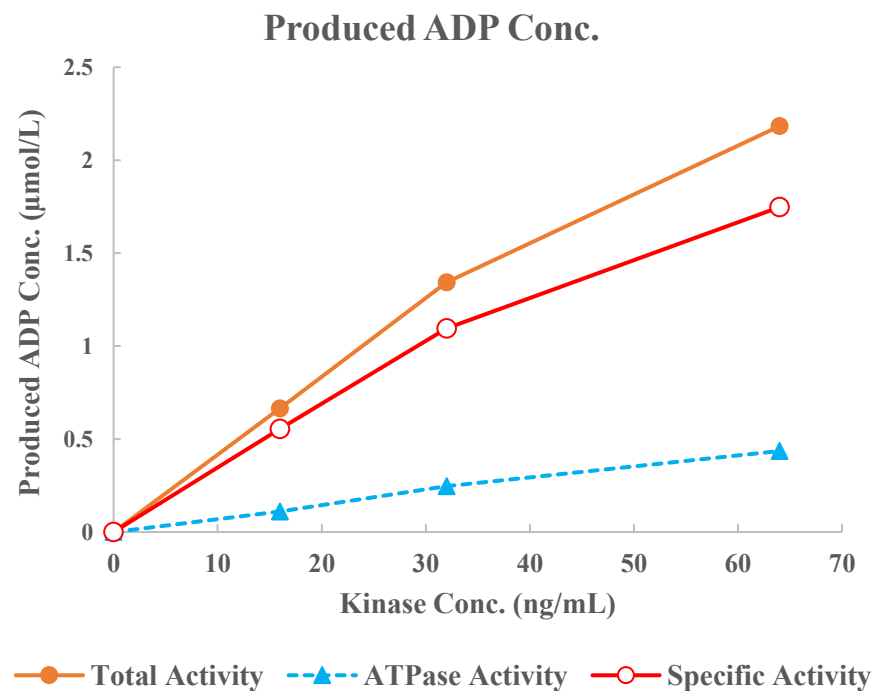
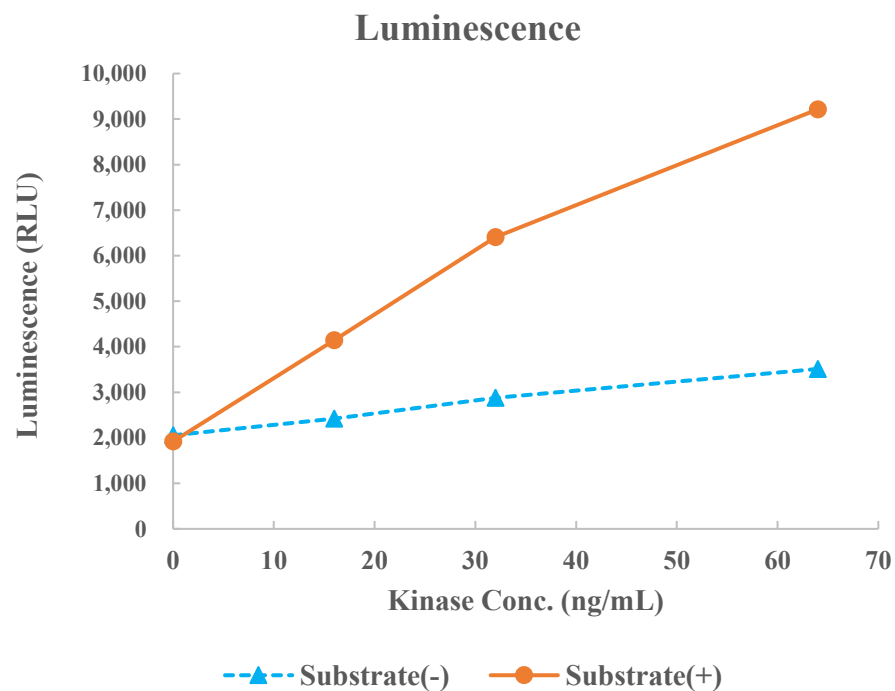


## CDK1(CDC2)/CycB1



|                 |                                    |
|-----------------|------------------------------------|
| Kinase          | CDK1(CDC2)/CycB1, 04-102           |
| Substrate       | Biotin-Modified Histone H1, BTSU13 |
| Substrate Conc. | 10 μmol/L                          |
| ATP Conc.       | 50 μmol/L                          |
| Metal           | Mg                                 |
| Metal Conc.     | 5 mmol/L                           |
| Additives       | -                                  |
| Reaction Time   | 1 hour                             |