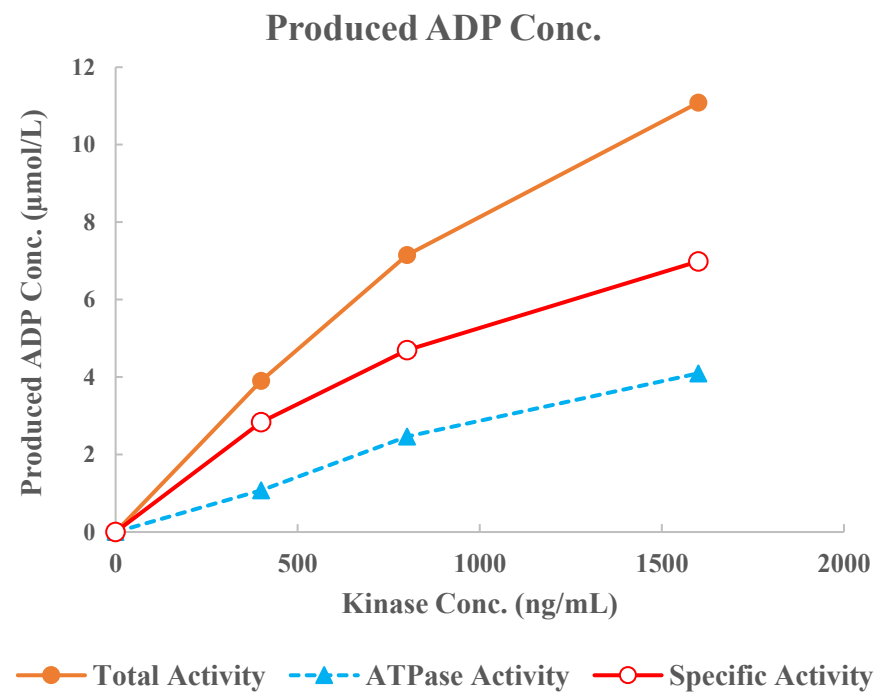
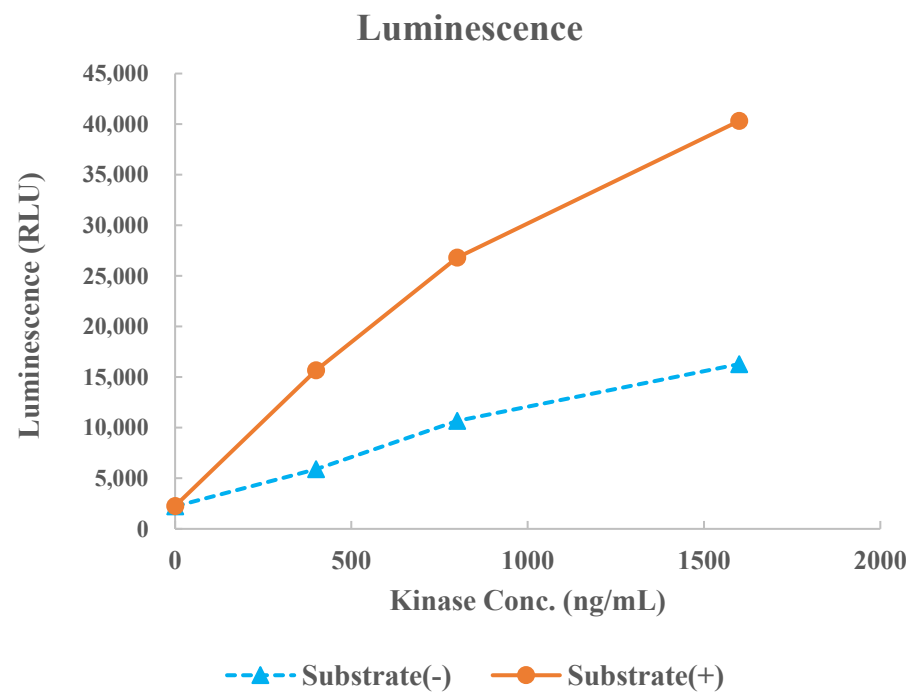


## NEK1



|                 |                             |
|-----------------|-----------------------------|
| Kinase          | NEK1, 05-123                |
| Substrate       | Biotin-CDK7 peptide, BTSU11 |
| Substrate Conc. | 10 μmol/L                   |
| ATP Conc.       | 75 μmol/L                   |
| Metal           | Mg                          |
| Metal Conc.     | 5 mmol/L                    |
| Additives       | -                           |
| Reaction Time   | 1 hour                      |