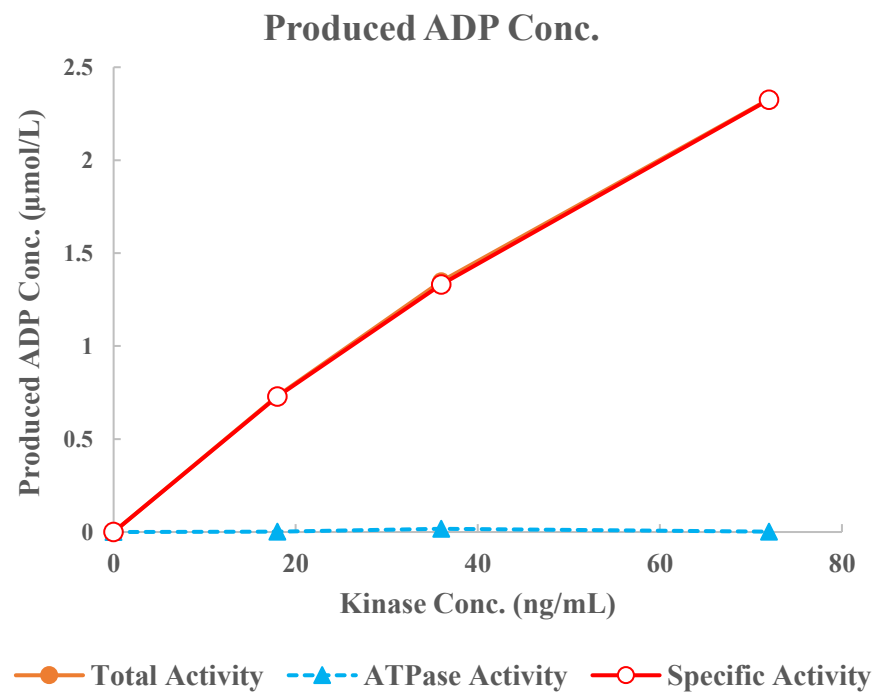
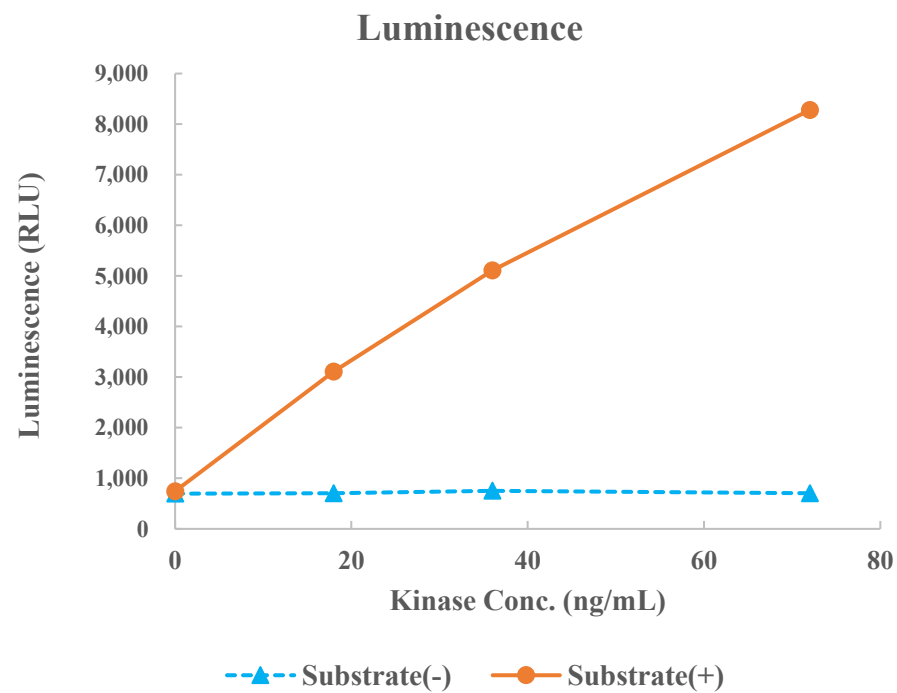


## EPHA1



|                 |                            |
|-----------------|----------------------------|
| Kinase          | EPHA1, 08-119              |
| Substrate       | Biotin-Blk/Lyntide, BTSU04 |
| Substrate Conc. | 10 μmol/L                  |
| ATP Conc.       | 25 μmol/L                  |
| Metal           | Mg                         |
| Metal Conc.     | 5 mmol/L                   |
| Additives       | -                          |
| Reaction Time   | 1 hour                     |