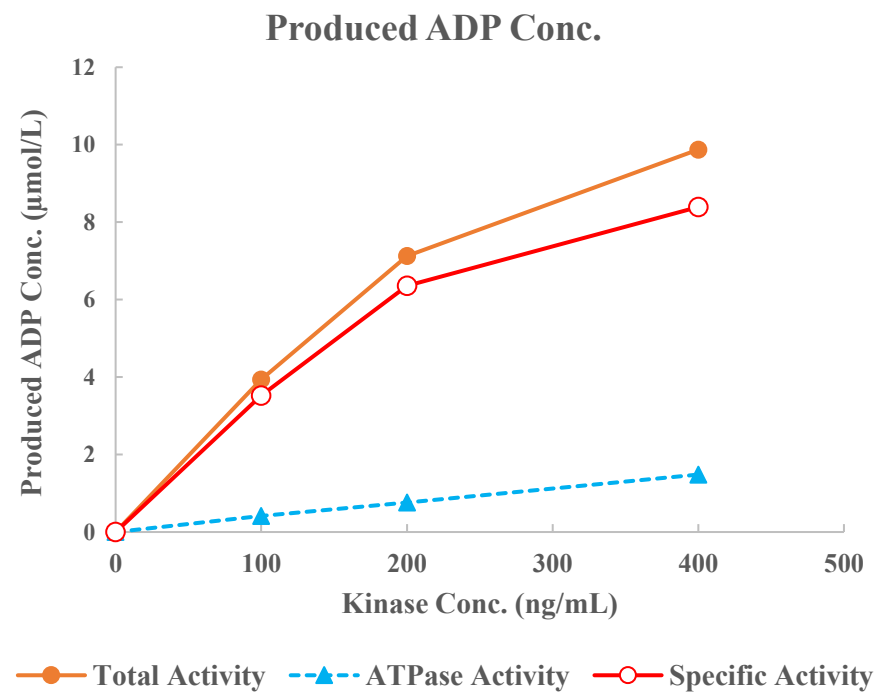
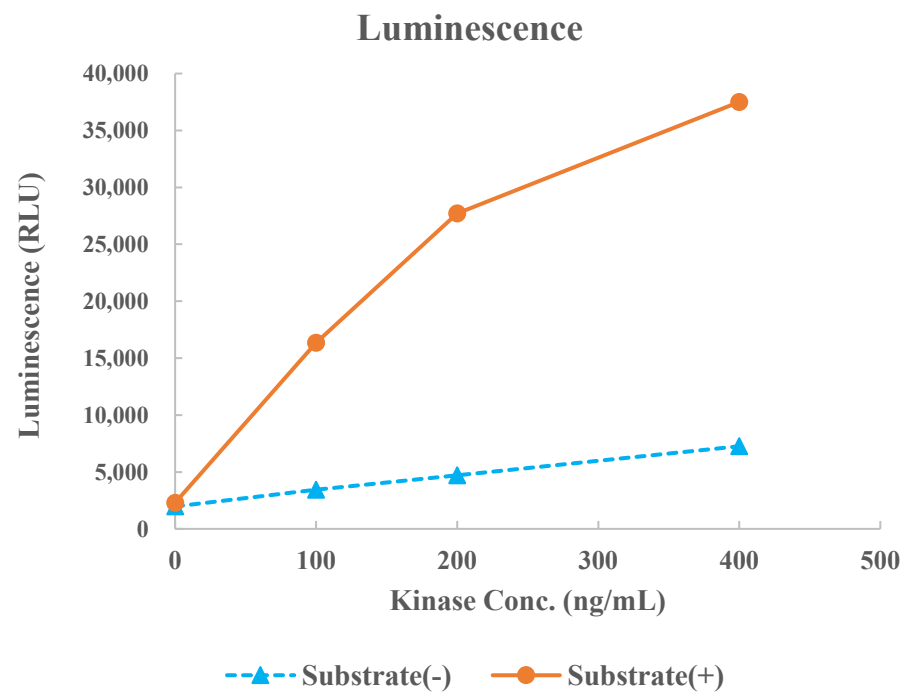


## TRKC(NTRK3)



|                 |                         |
|-----------------|-------------------------|
| Kinase          | TRKC(NTRK3), 08-197     |
| Substrate       | Biotin-Srcptide, BTSU01 |
| Substrate Conc. | 10 μmol/L               |
| ATP Conc.       | 50 μmol/L               |
| Metal           | Mg                      |
| Metal Conc.     | 5 mmol/L                |
| Additives       | -                       |
| Reaction Time   | 1 hour                  |