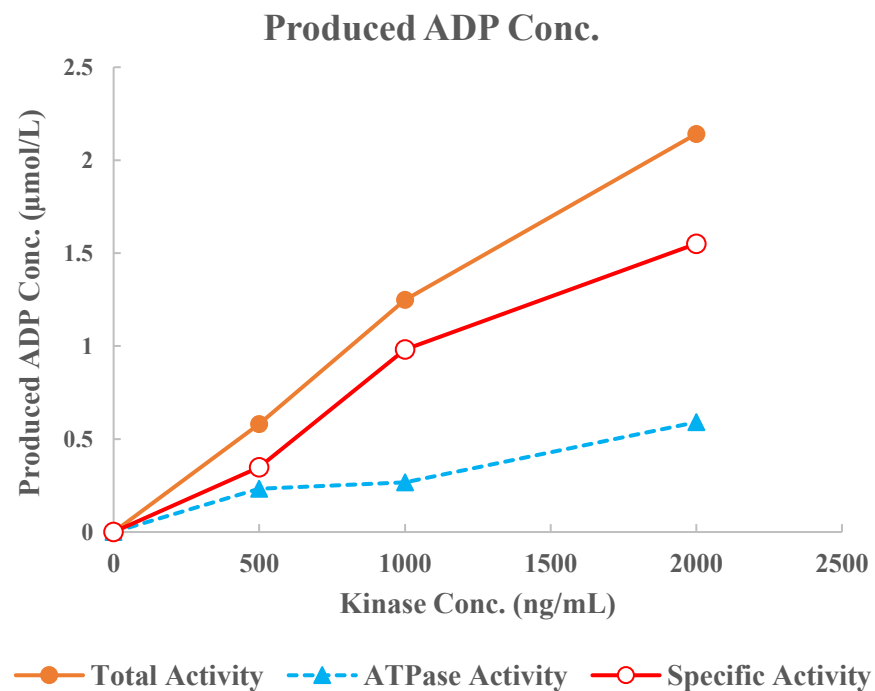
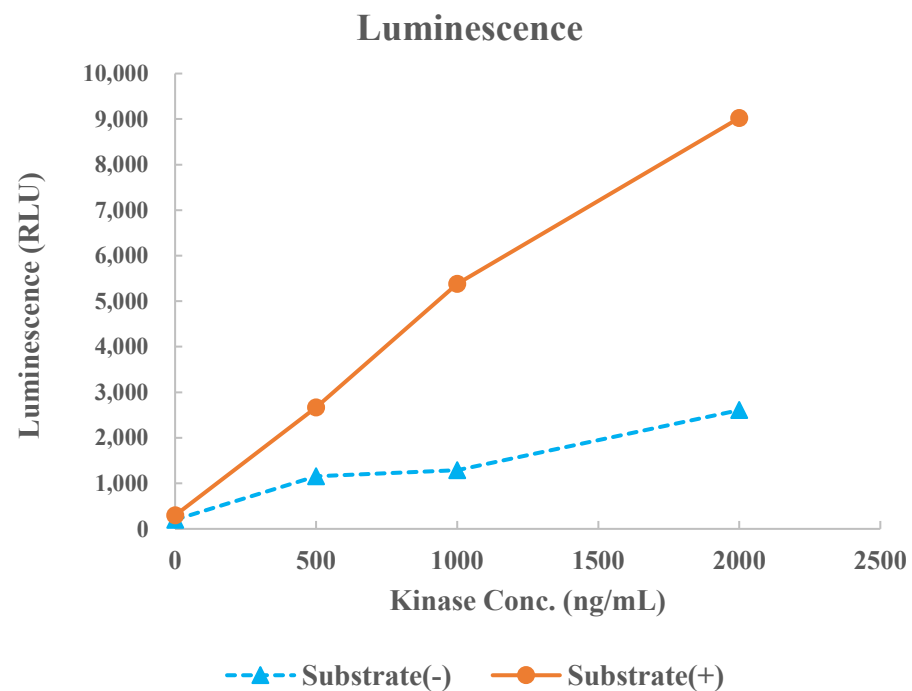


## MRCK $\beta$ (CDC42BPB)



|                 |                                 |
|-----------------|---------------------------------|
| Kinase          | MRCK $\beta$ (CDC42BPB), 01-108 |
| Substrate       | Biotin-DAPK1tide, BTSU17        |
| Substrate Conc. | 10 $\mu$ mol/L                  |
| ATP Conc.       | 10 $\mu$ mol/L                  |
| Metal           | Mg                              |
| Metal Conc.     | 5 mmol/L                        |
| Additives       | -                               |
| Reaction Time   | 1 hour                          |