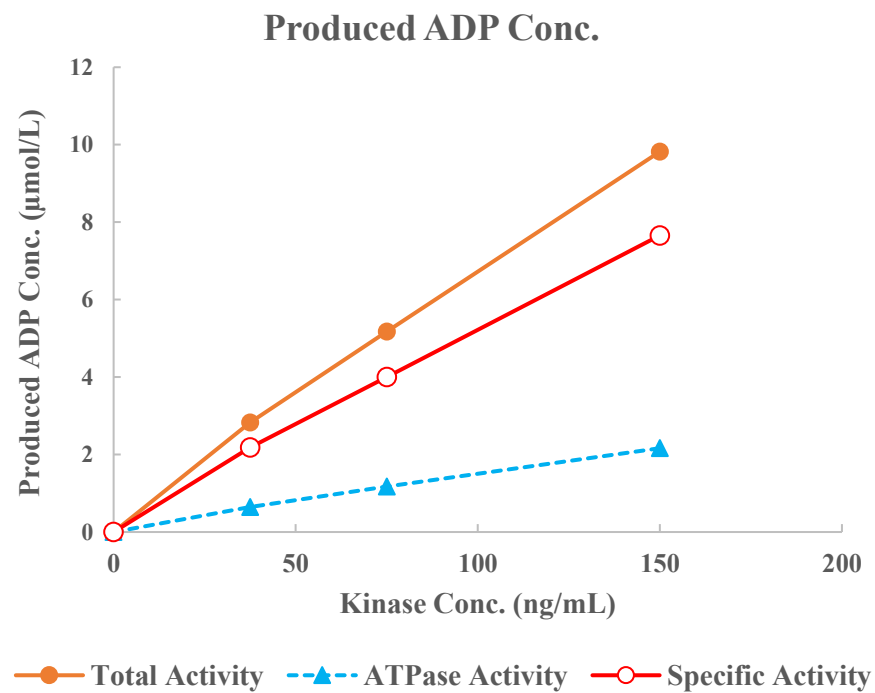
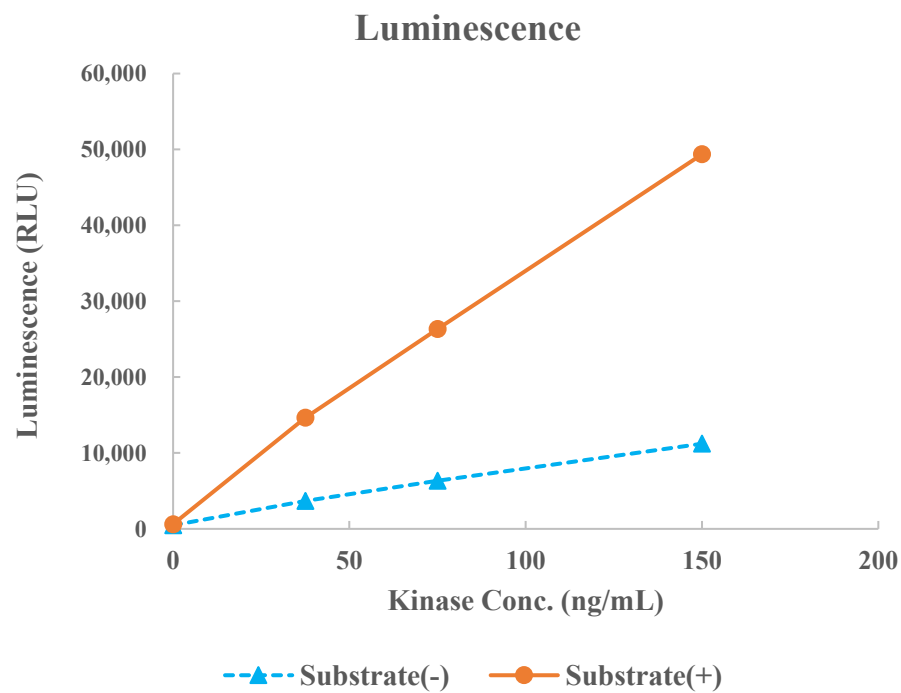


## p38δ(MAPK13)



|                 |                                 |
|-----------------|---------------------------------|
| Kinase          | p38δ(MAPK13), 04-154            |
| Substrate       | Biotin-Modified Erktide, BTSU03 |
| Substrate Conc. | 50 μmol/L                       |
| ATP Conc.       | 10 μmol/L                       |
| Metal           | Mg                              |
| Metal Conc.     | 5 mmol/L                        |
| Additives       | -                               |
| Reaction Time   | 1 hour                          |