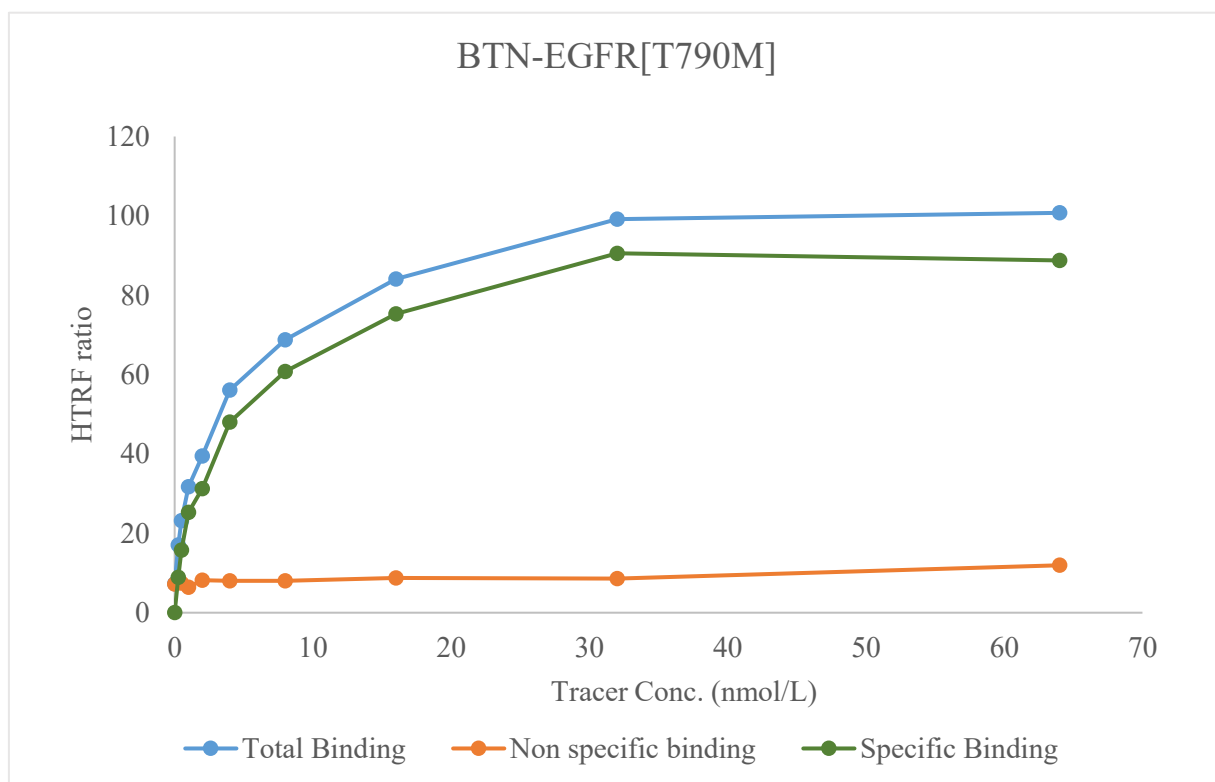


Kinase **BTN-EGFR[T790M]**

Cat.# 08-494-20N



|                         |                                                  |
|-------------------------|--------------------------------------------------|
| Tracer                  | revvity Inc.<br>Staurosporine-Red, (#62KB01REDC) |
| KD                      | 4.2 nmol/L                                       |
| Kinase Conc.            | 5 nmol/L                                         |
| Eu labeled-streptavidin | 5 nmol/L                                         |
| Metal                   | Mg, Mn                                           |
| Metal Conc.             | 5, 1 mmol/L                                      |
| Additives               | -                                                |
| Reaction Time           | 1 Hour                                           |