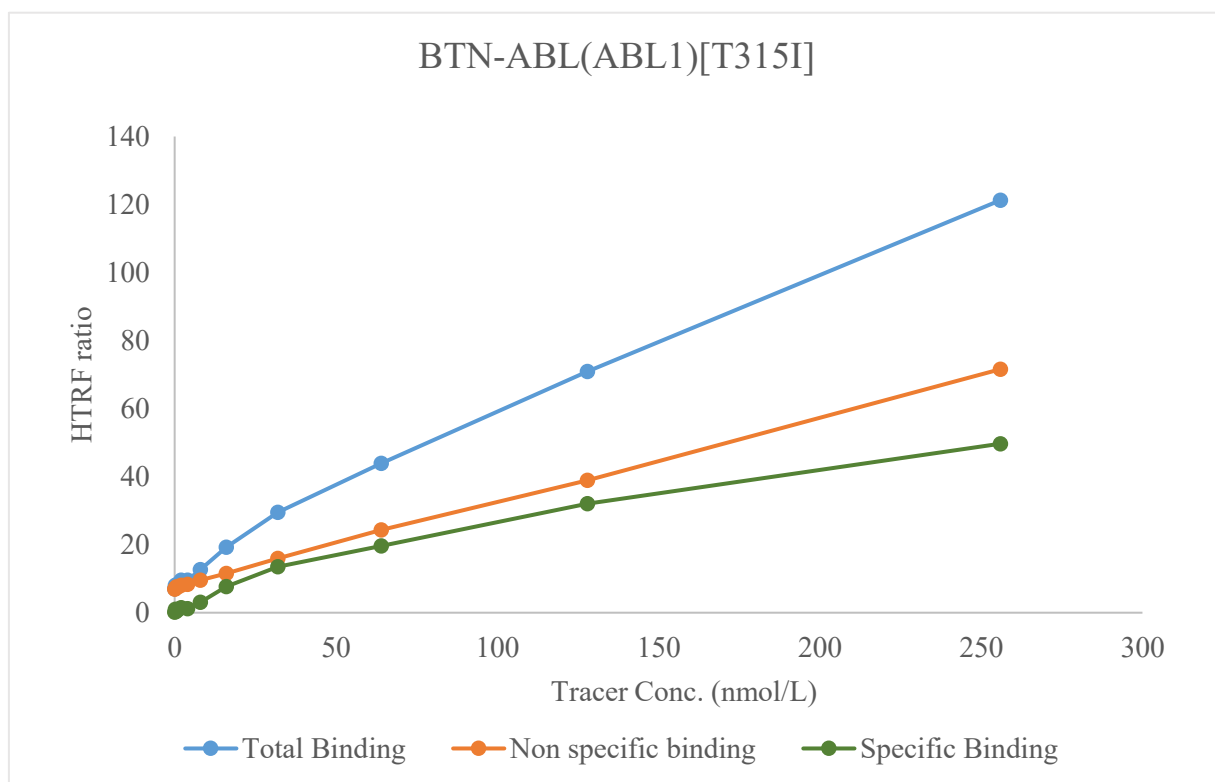


Kinase **BTN-ABL(ABL1)[T315I]**

Cat.# 08-493-20N



|                         |                                              |
|-------------------------|----------------------------------------------|
| Tracer                  | revvity Inc.<br>Dasatinib-Red, (#62KB02REDC) |
| KD                      | 210 nmol/L                                   |
| Kinase Conc.            | 5 nmol/L                                     |
| Eu labeled-streptavidin | 5 nmol/L                                     |
| Metal                   | Mg                                           |
| Metal Conc.             | 5 mmol/L                                     |
| Additives               | -                                            |
| Reaction Time           | 1 Hour                                       |