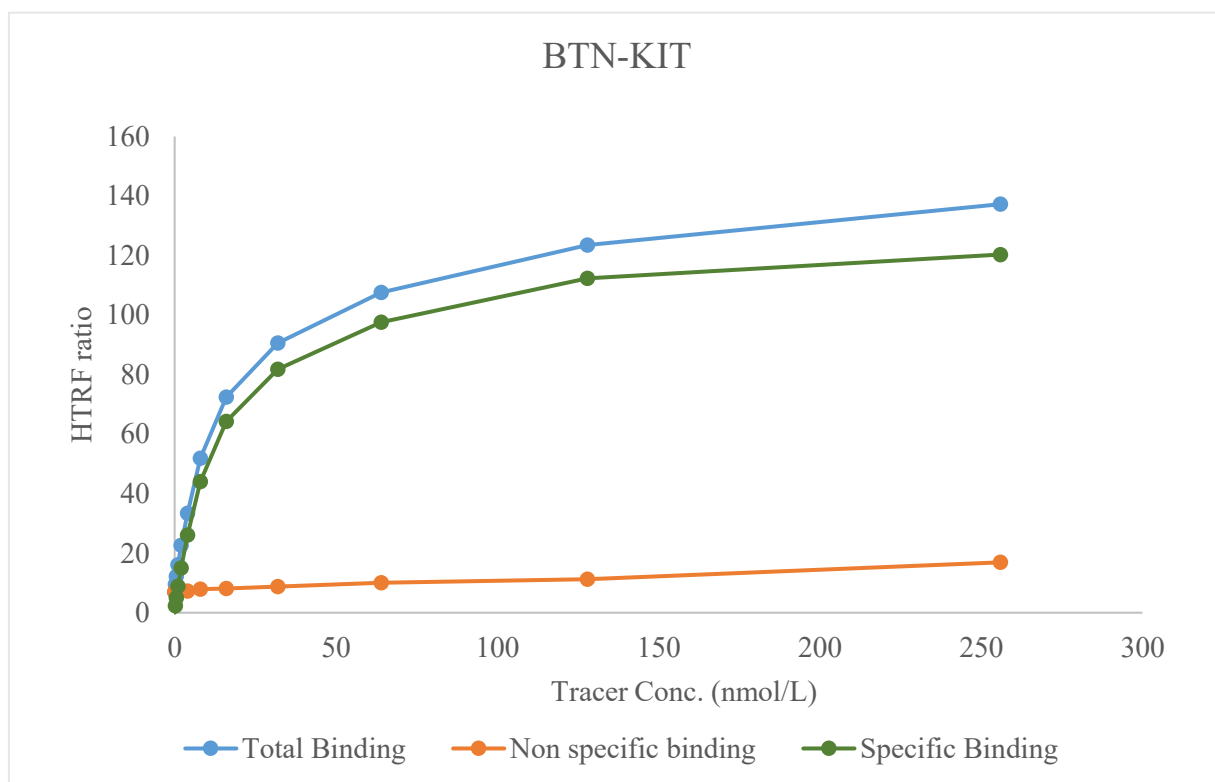


Kinase **BTN-KIT**

Cat.# 08-456-21N



|                         |                                              |
|-------------------------|----------------------------------------------|
| Tracer                  | revvity Inc.<br>Sunitinib-Red, (#62KB03REDC) |
| KD                      | 15 nmol/L                                    |
| Kinase Conc.            | 5 nmol/L                                     |
| Eu labeled-streptavidin | 5 nmol/L                                     |
| Metal                   | Mg                                           |
| Metal Conc.             | 5 mmol/L                                     |
| Additives               | 0.025 mmol/L Sodium orthovanadate            |
| Reaction Time           | 1 Hour                                       |