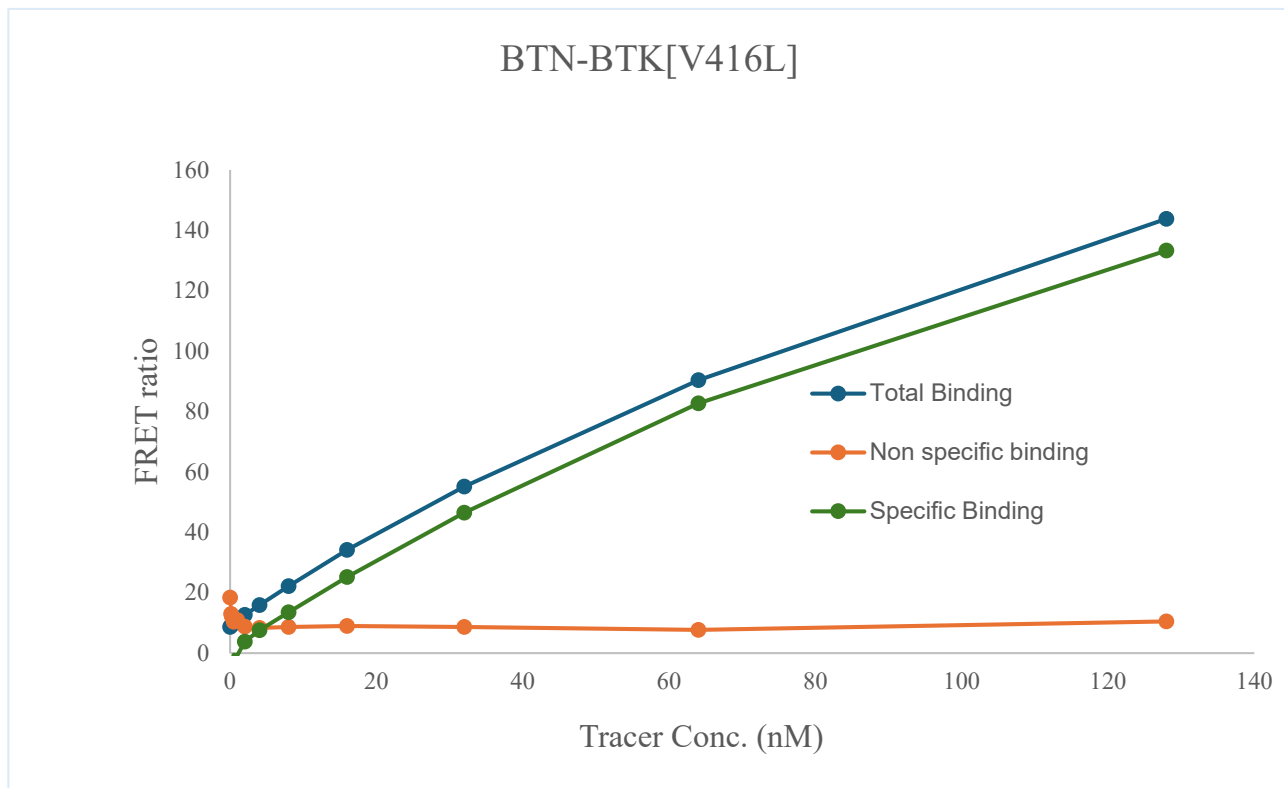


Kinase **BTN-BTK[V416L]**  
 Cat.# 08-706-20N  
 Tracer Revvity Inc.  
 Staurosporine-Red (#62KB02REDC)



|                         |     |   |
|-------------------------|-----|---|
| KD (nM)                 | 217 |   |
| Kinase conc.            |     | 5 |
| Eu labeled-streptavidin |     | 5 |

Follow the guide to set up your HTRF® binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF® kinase binding

