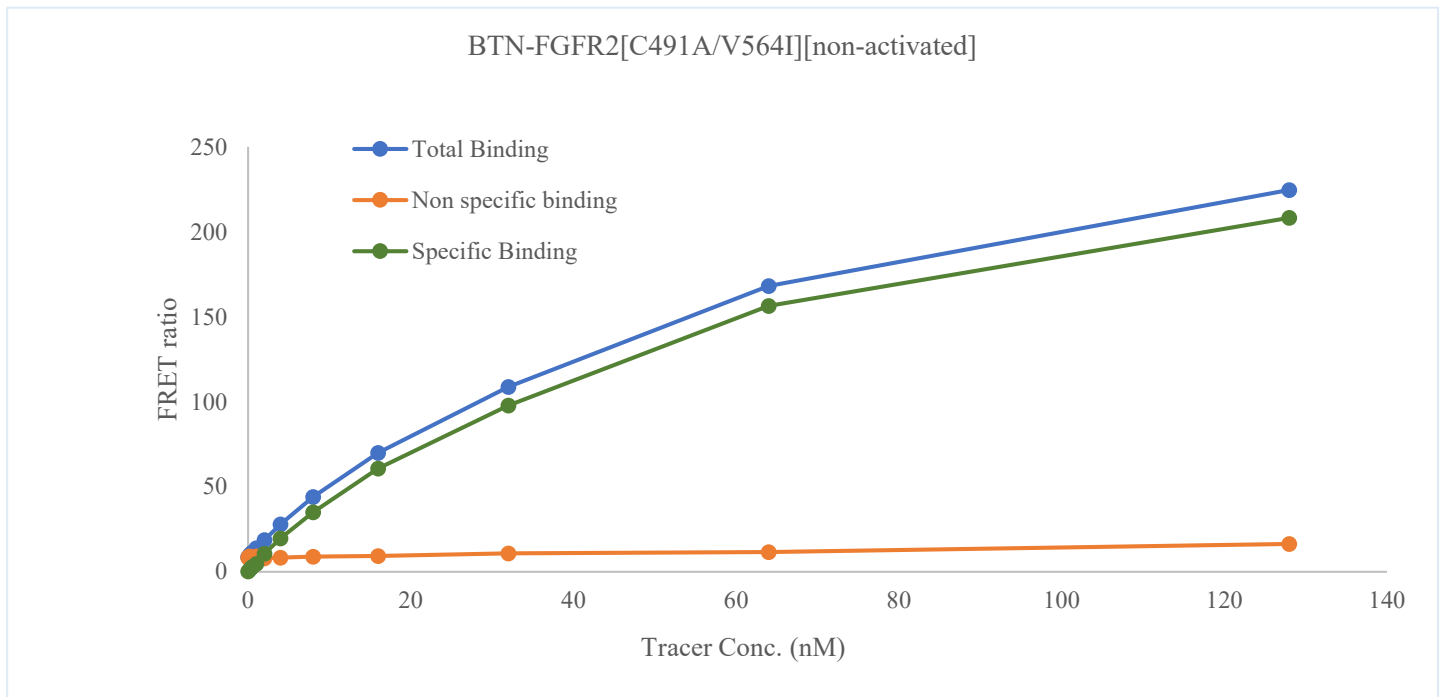


Kinase **BTN-FGFR2[C491A/V564I][non-activated]**
 Cat.# 08-905-23N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB01REDC)



KD (nM)	70.2
Kinase conc.	5nM
Eu labeled-streptavidin	5nM

Follow the guide to set up your HTRF® binding assay



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

