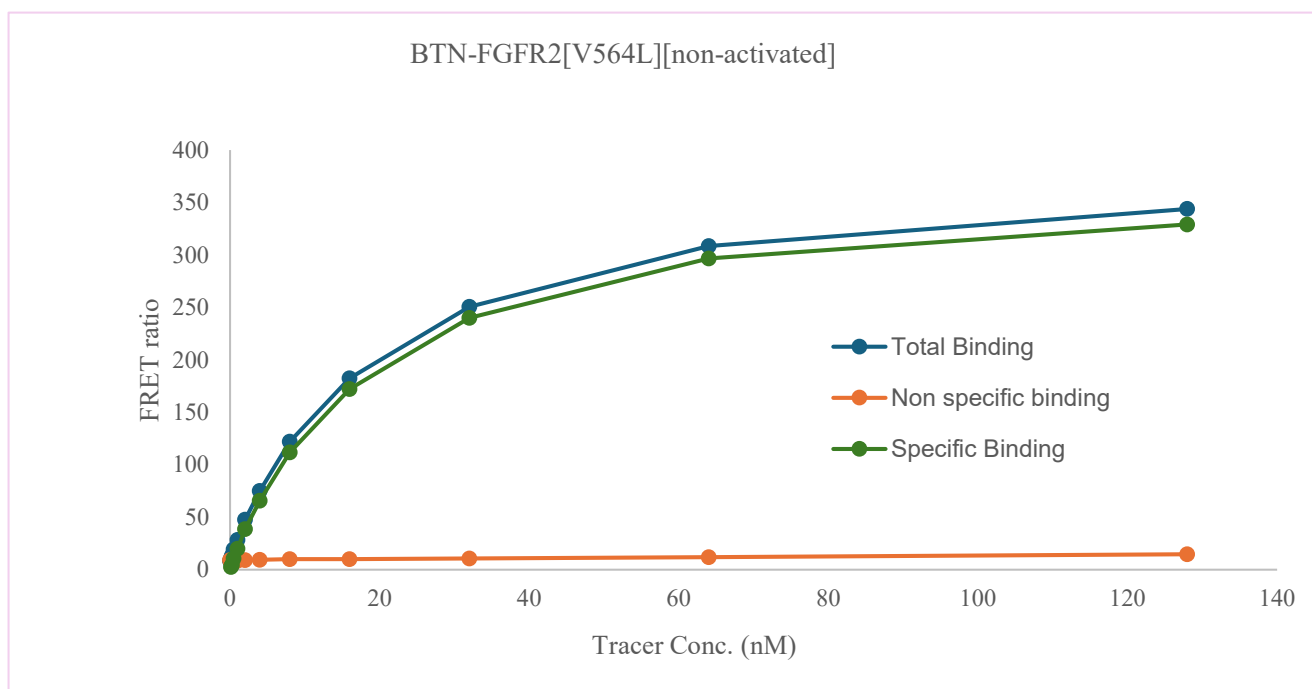


Kinase **BTN-FGFR2[V564L][non-activated]**
 Cat.# 08-904-23N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB02REDC)



KD (nM)	18.9
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

