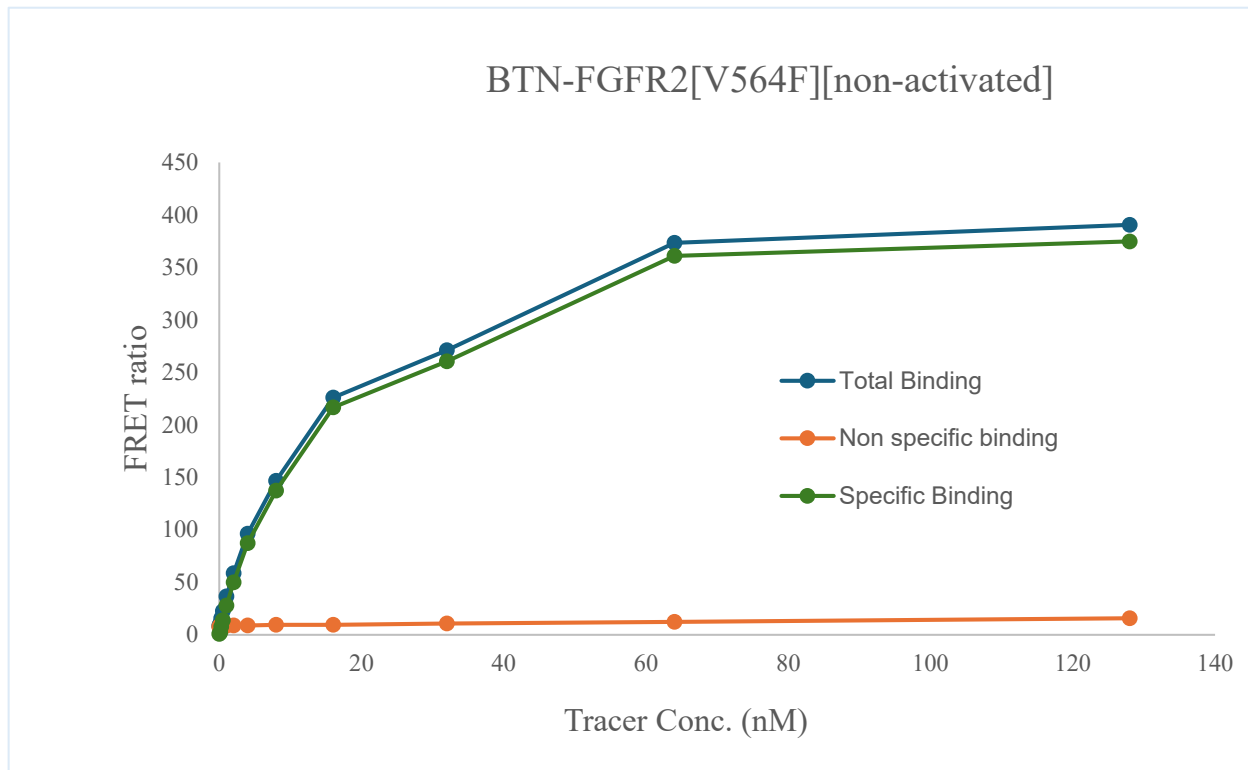


Kinase **BTN-FGFR2[V564F][non-activated]**
 Cat.# 08-910-23N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB01REDC)



KD (nM) 16.9
 Kinase conc. 5 nM
 Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF® binding assay

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

