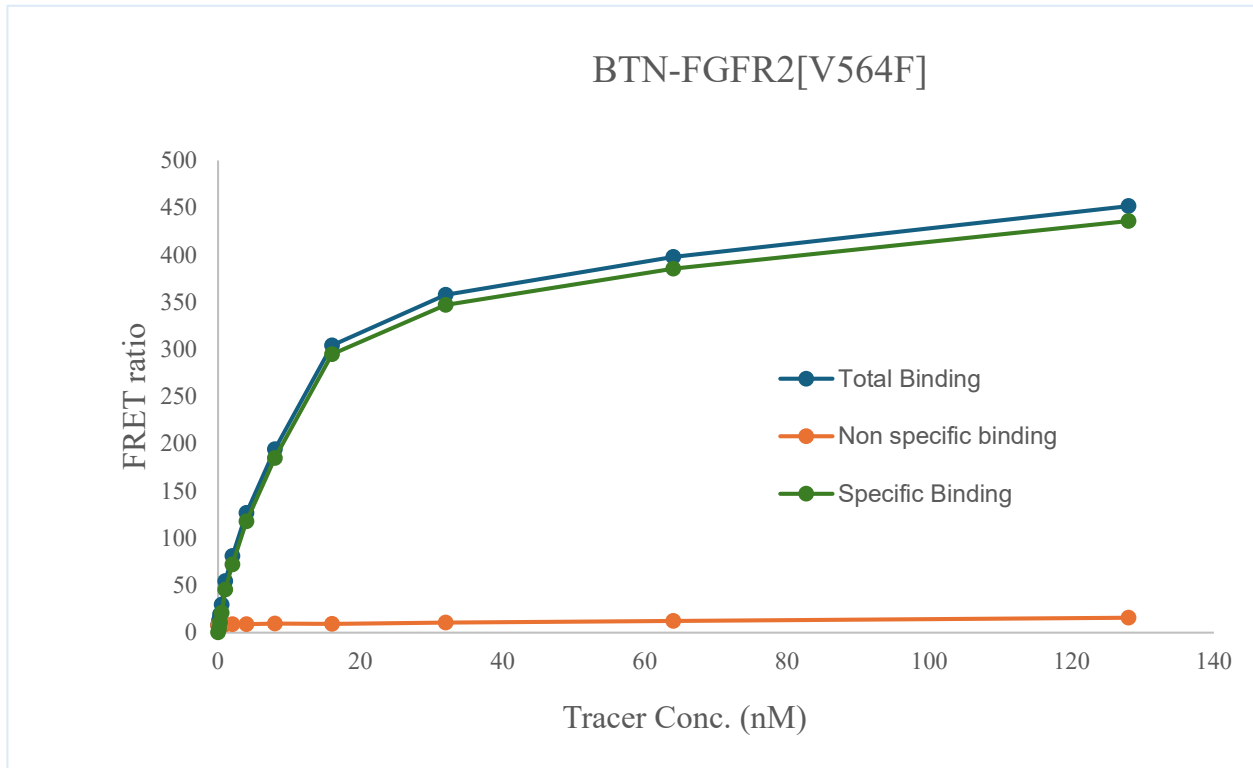


Kinase **BTN-FGFR2[V564F]**  
 Cat.# 08-910-20N  
 Tracer Revvity Inc.  
 Staurosporine-Red (#62KB01REDC)



KD (nM) 11  
 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

