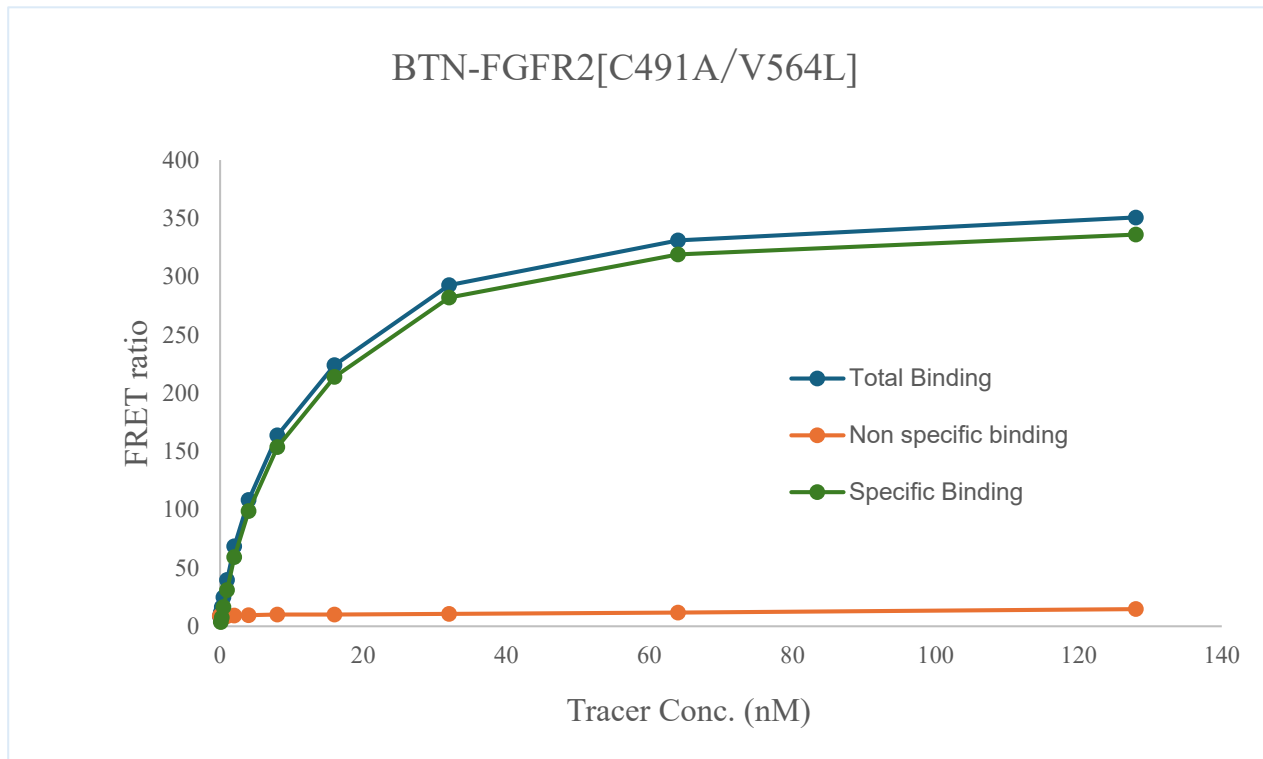


Kinase **BTN-FGFR2[C491A/V564L]**  
 Cat.# 08-906-20N  
 Tracer Revvity Inc.  
 Staurosporine-Red (#62KB02REDC)



KD (nM) 11.1  
 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

