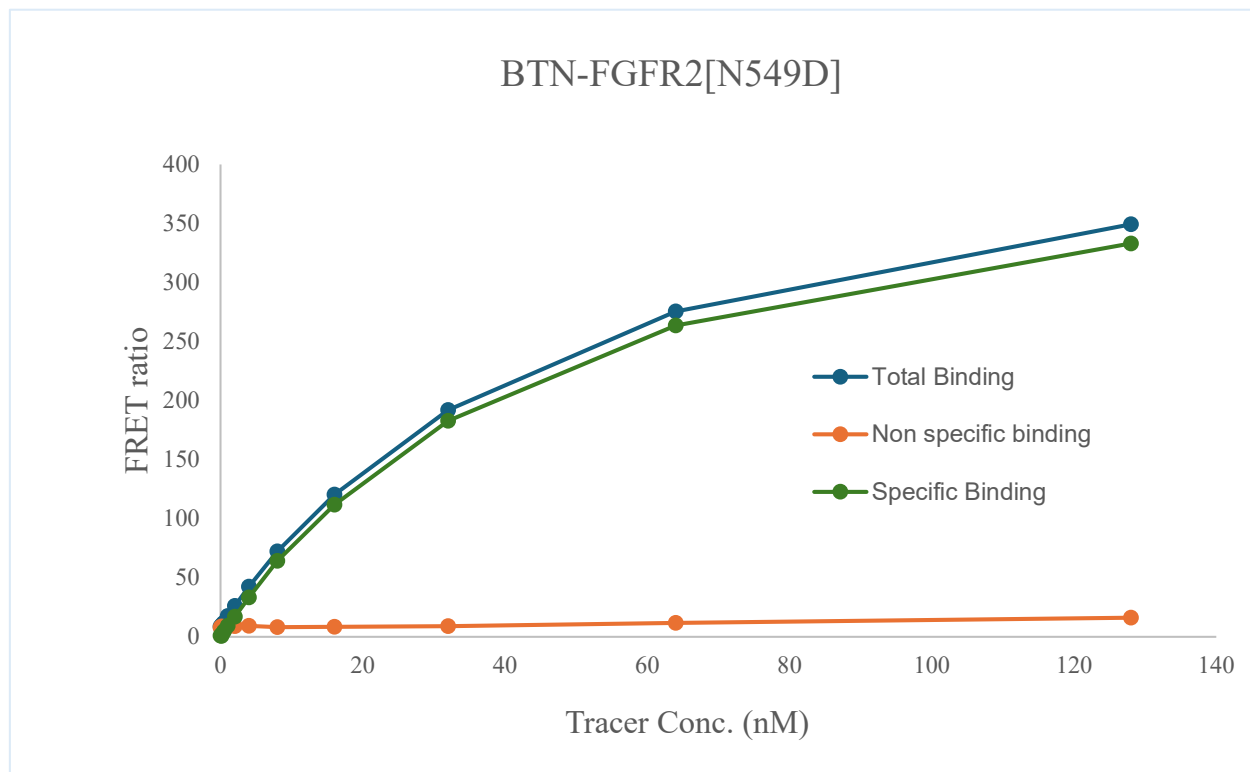


Kinase **BTN-FGFR2[N549D]**
 Cat.# 08-908-20N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB02REDC)



KD (nM) 50

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

