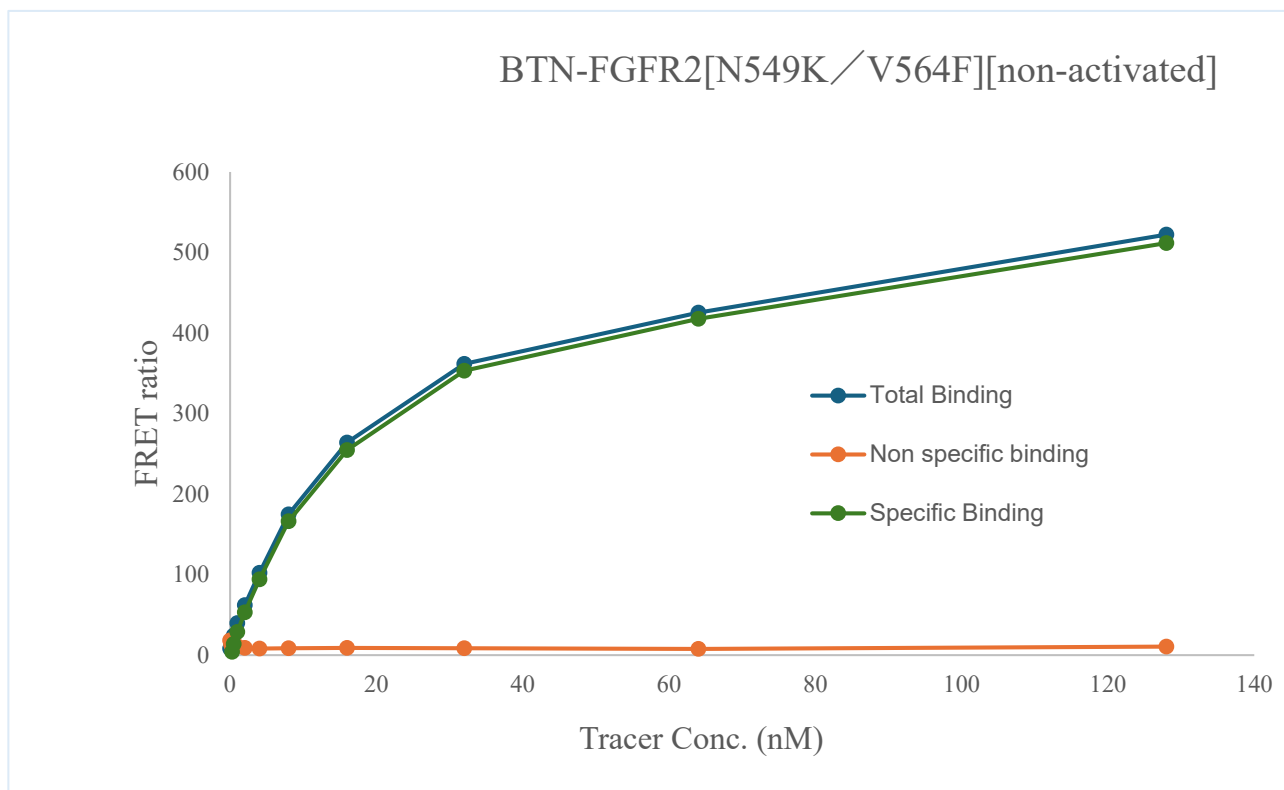


Kinase **BTN-FGFR2[N549K/V564F][non-activated]**
 Cat.# 08-925-23N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB02REDC)



KD (nM)	19.9	
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

