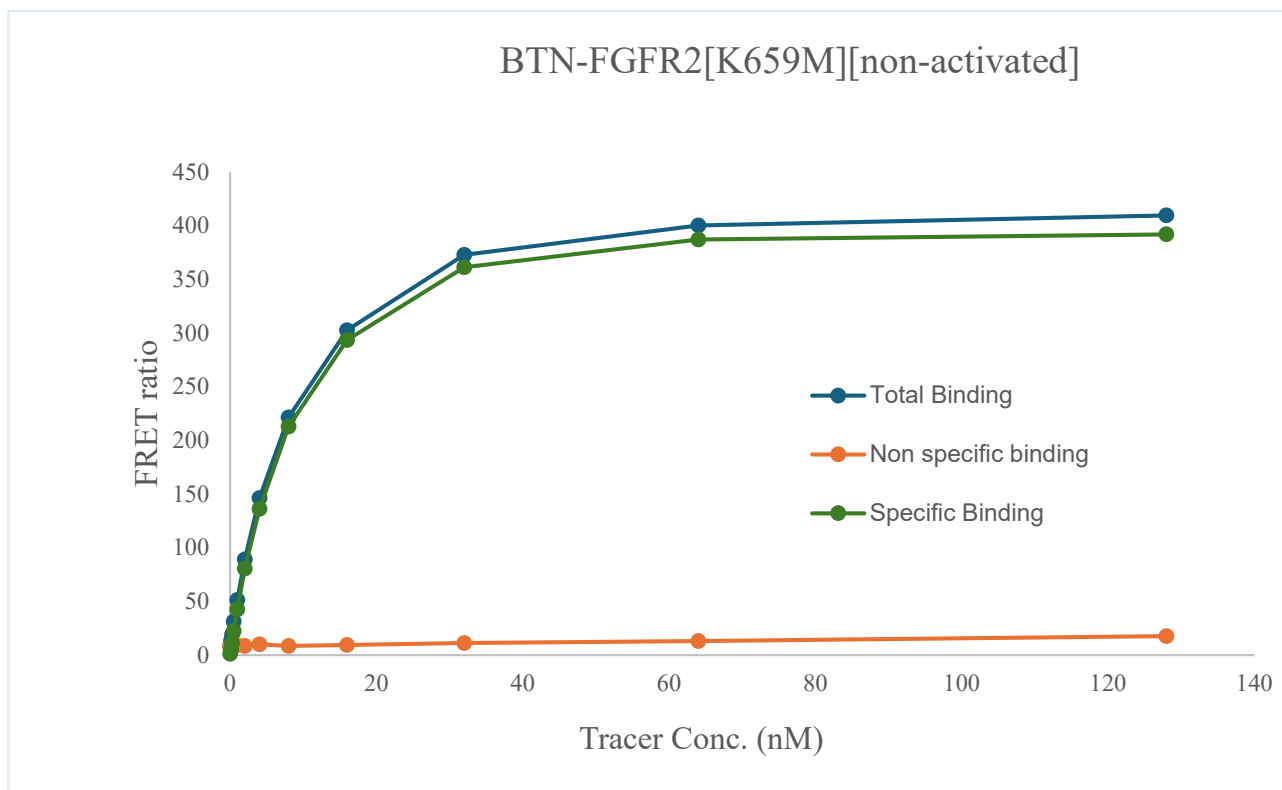


Kinase **BTN-FGFR2[K659M][non-activated]**

Cat.# 08-927-23N

Tracer Revvity Inc.

Staurosporine-Red (#62KB02REDC)



KD (nM) 8.3

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

