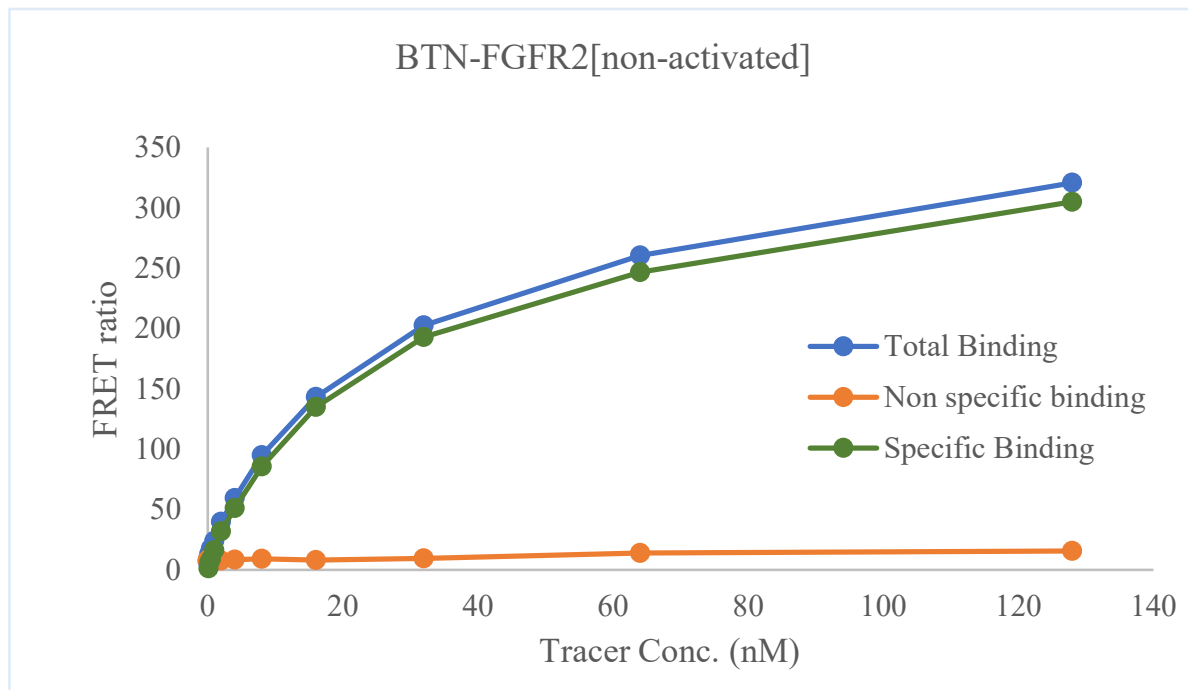


Kinase **BTN-FGFR2[non-activated]**

Cat.# 08-434-23N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM) 27.1

Kinase conc. 5nM

Eu labeled-streptavidin 5nM

Follow the guide to set up your HTRF® binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF® kinase binding

