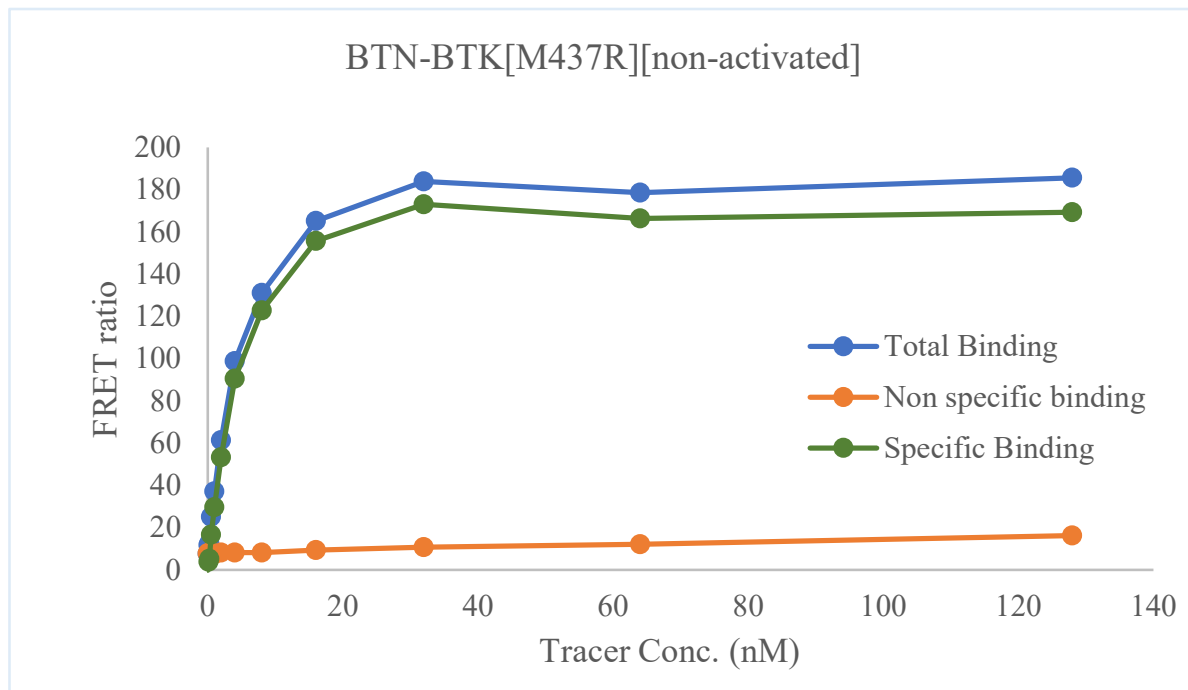


Kinase **BTN-BTK[M437R][non-activated]**

Cat.# 08-705-23N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM) 4.38

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

