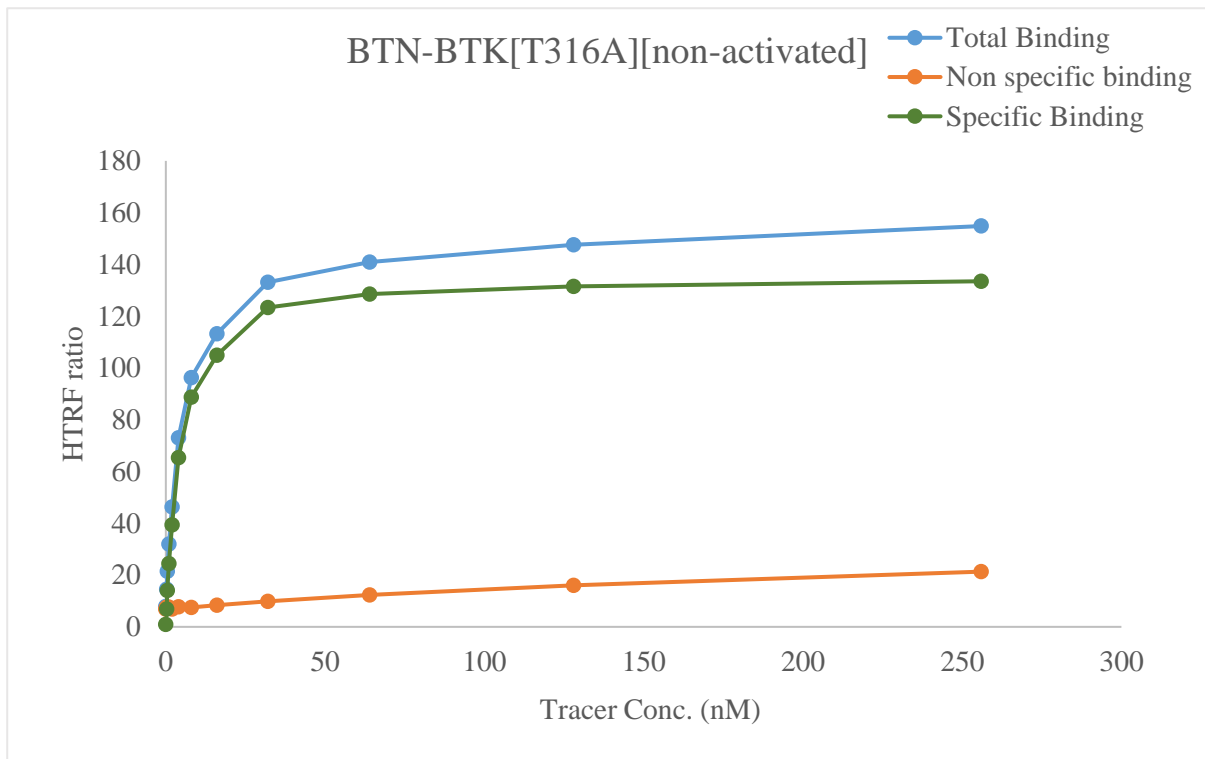


Kinase **BTN-BTK[T316A][non-activated]**

Cat.# 08-418-23N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 4.7

Kinase conc. 5 nM

Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF<sup>®</sup> binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF<sup>®</sup> kinase binding

