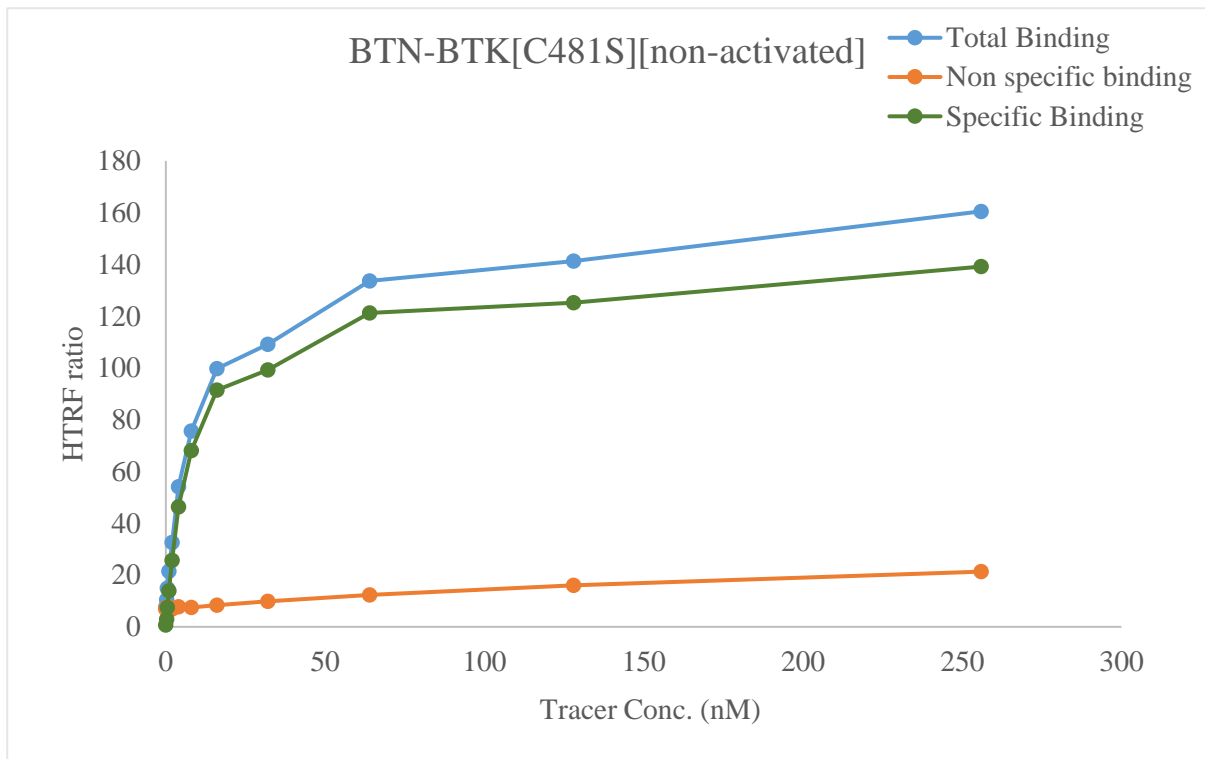


Kinase **BTN-BTK[C481S][non-activated]**  
 Cat.# 08-417-23N  
 Tracer revvity Inc.  
 Staurosporine-Red (#62KB01REDC)



KD (nM) 8.3

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

