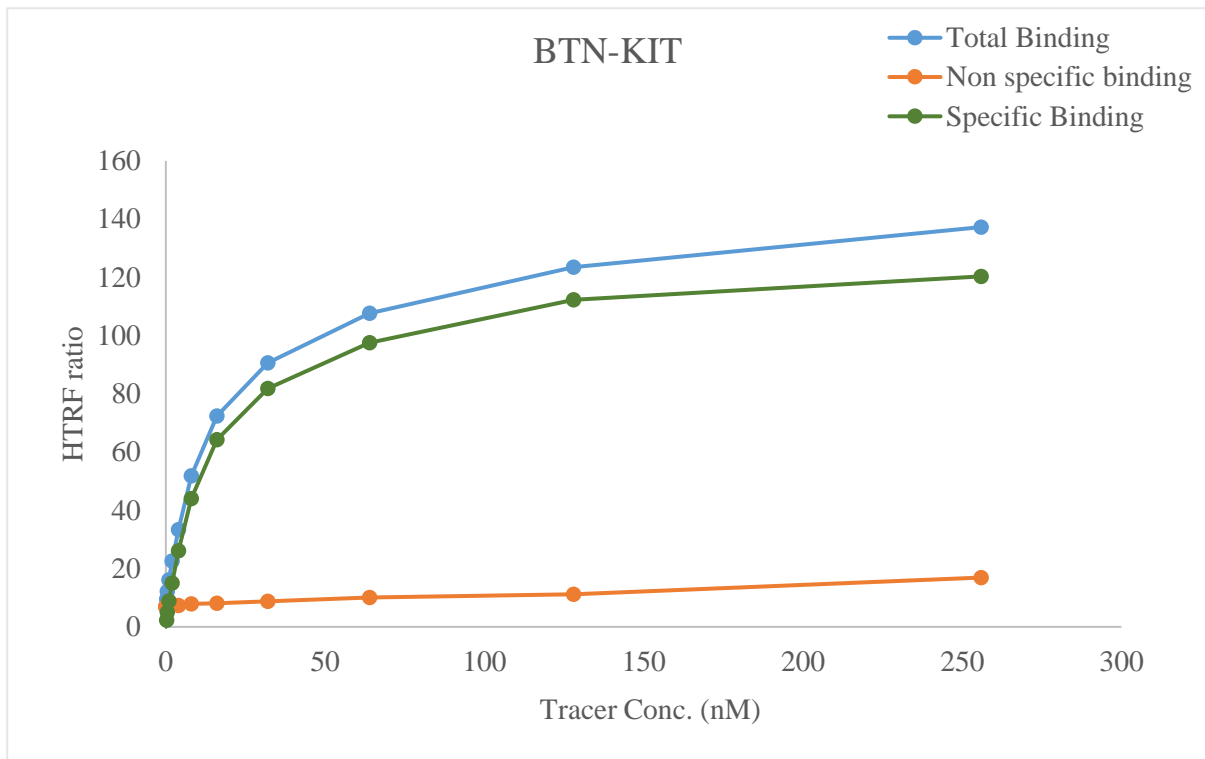


Kinase **BTN-KIT**  
 Cat.# 08-456-21N  
 Tracer revvity Inc.  
 Sunitinib-Red (#62KB03REDC)

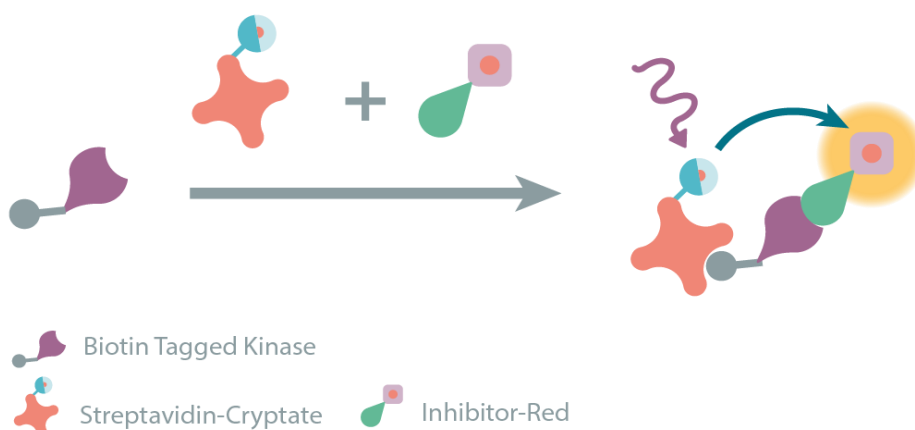


KD (nM) 15

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

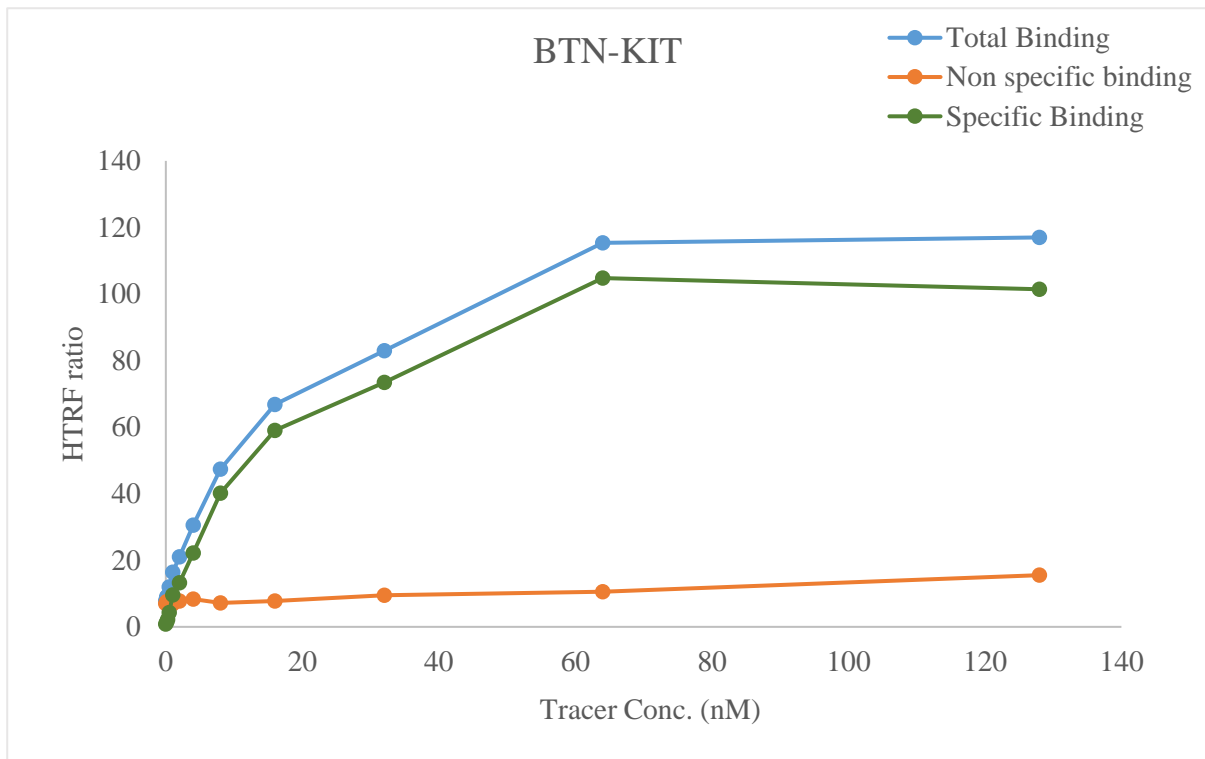


Kinase **BTN-KIT**

Cat.# 08-456-21N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 17

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

