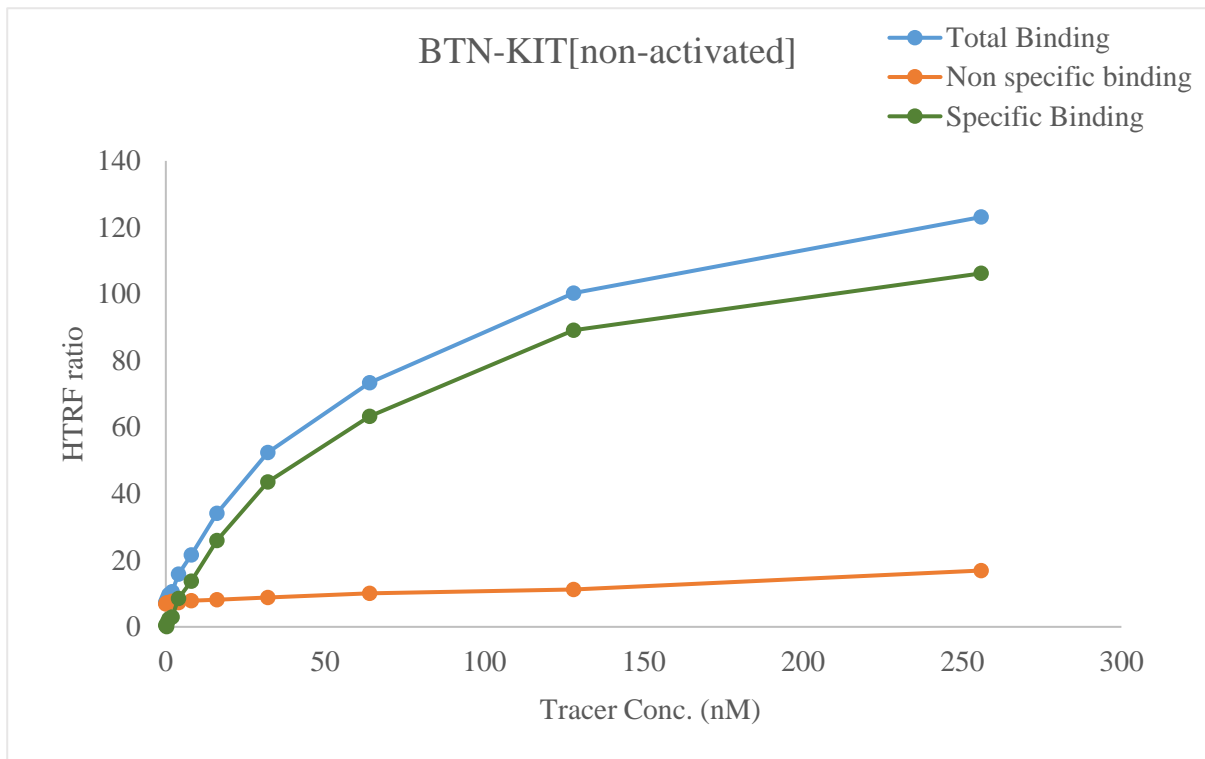


Kinase **BTN-KIT[non-activated]**
 Cat.# 08-456-23N
 Tracer revvity Inc.
 Sunitinib-Red (#62KB03REDC)

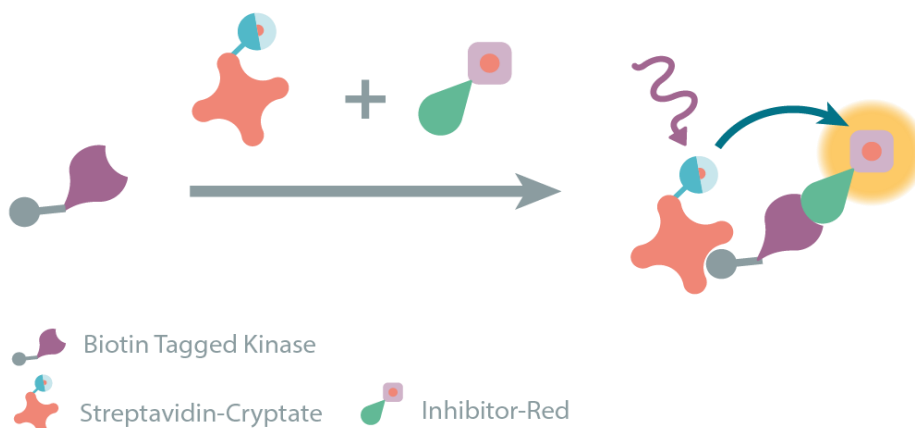


KD (nM) 68

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

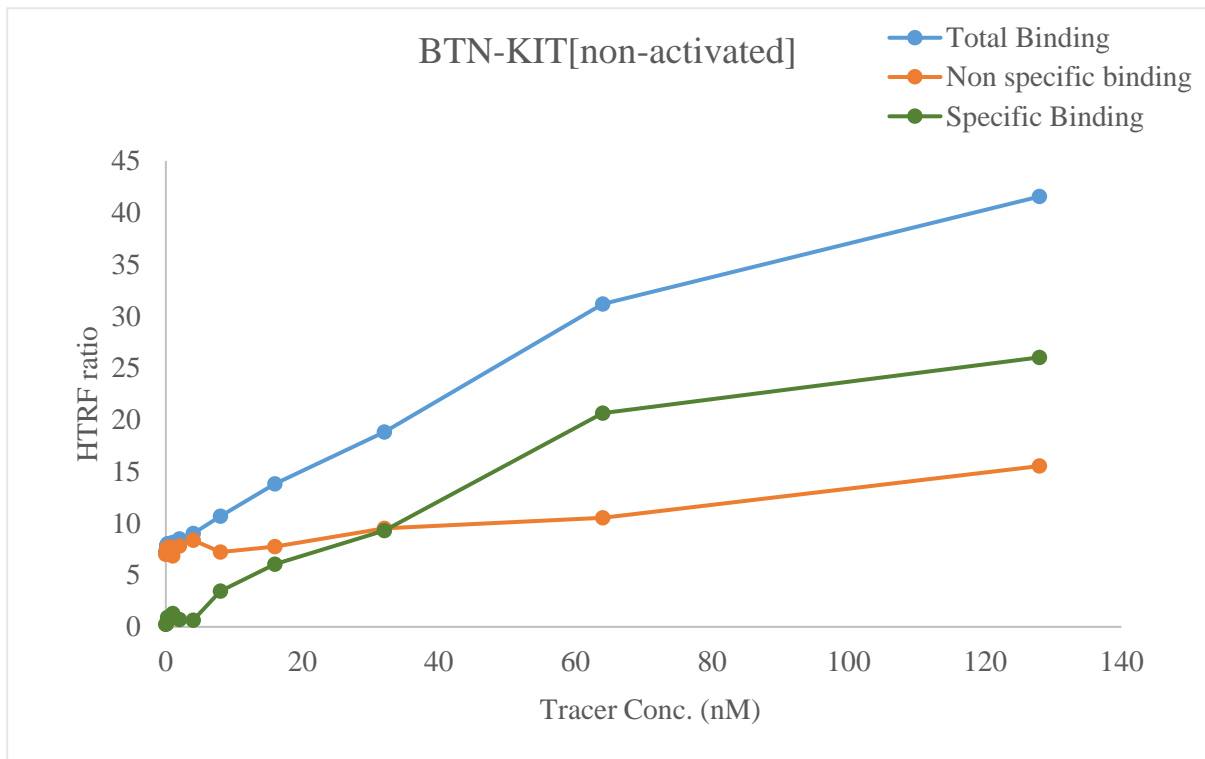


Kinase **BTN-KIT[non-activated]**

Cat.# 08-456-23N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 130

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

