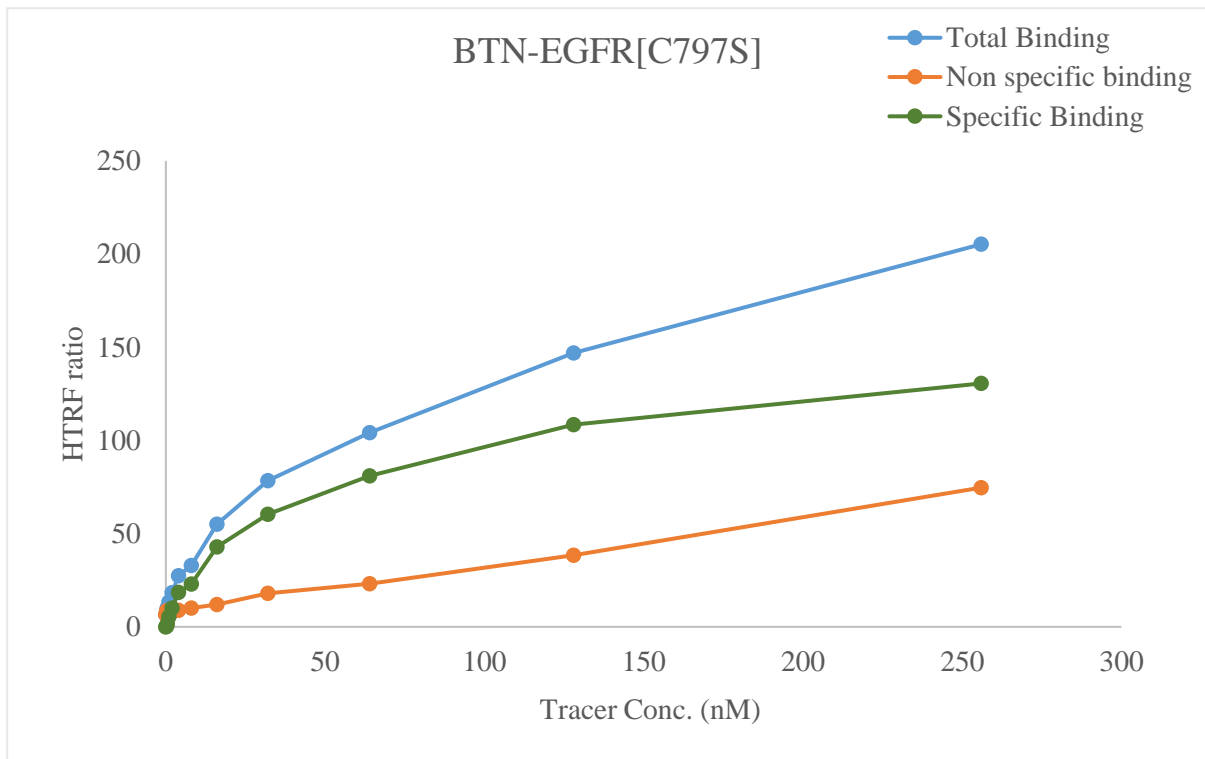


Kinase **BTN-EGFR[C797S]**
 Cat.# 08-761-20N
 Tracer revvity Inc.
 Dasatinib-Red (#62KB02REDC)

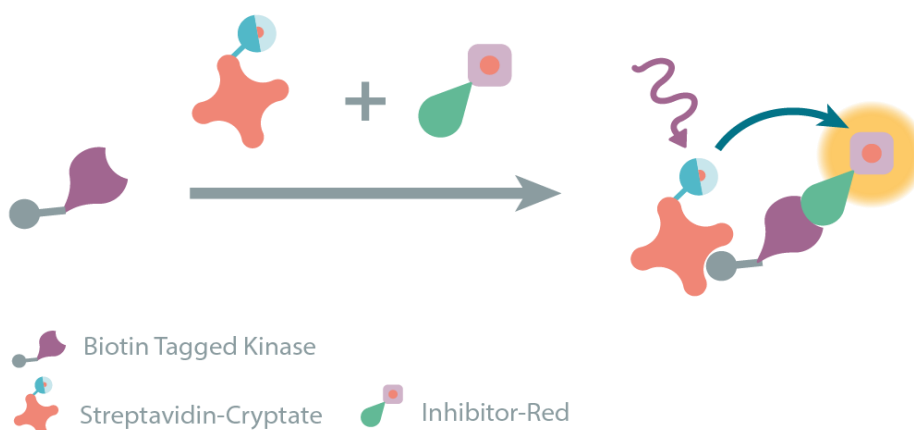


KD (nM) 46

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

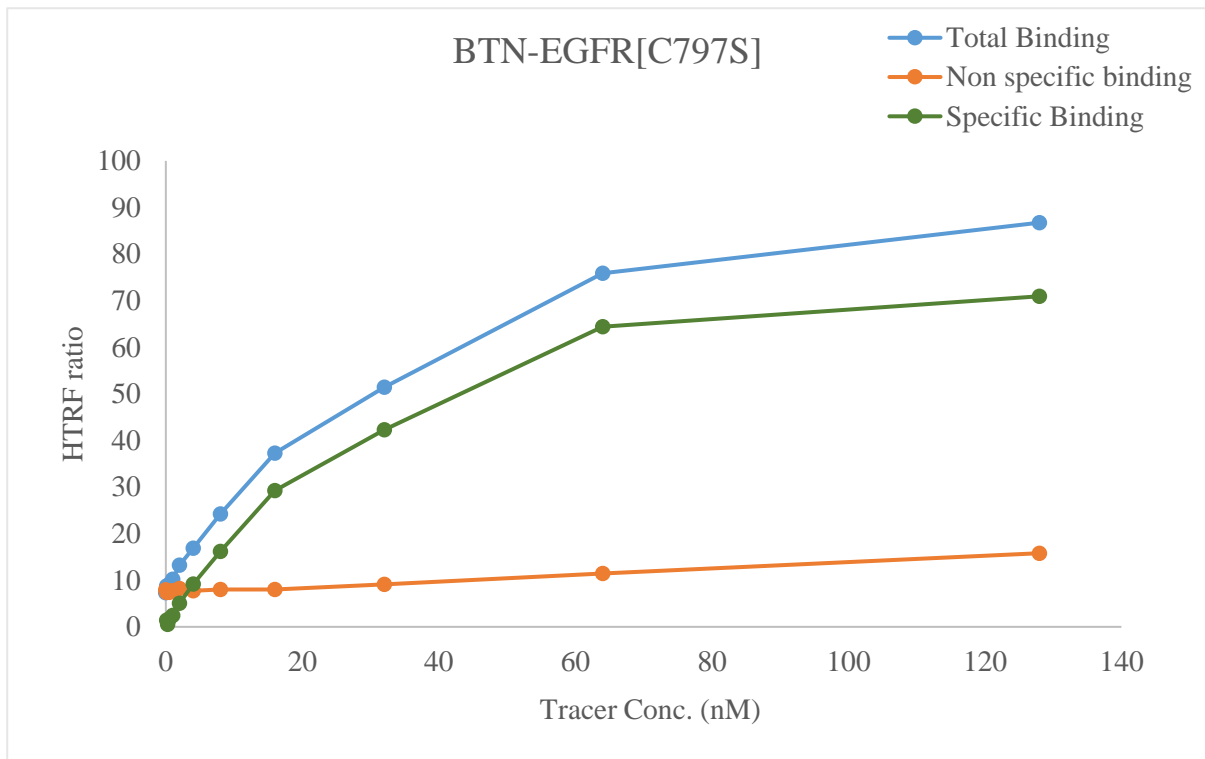


Kinase **BTN-EGFR[C797S]**

Cat.# 08-761-20N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 37

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

