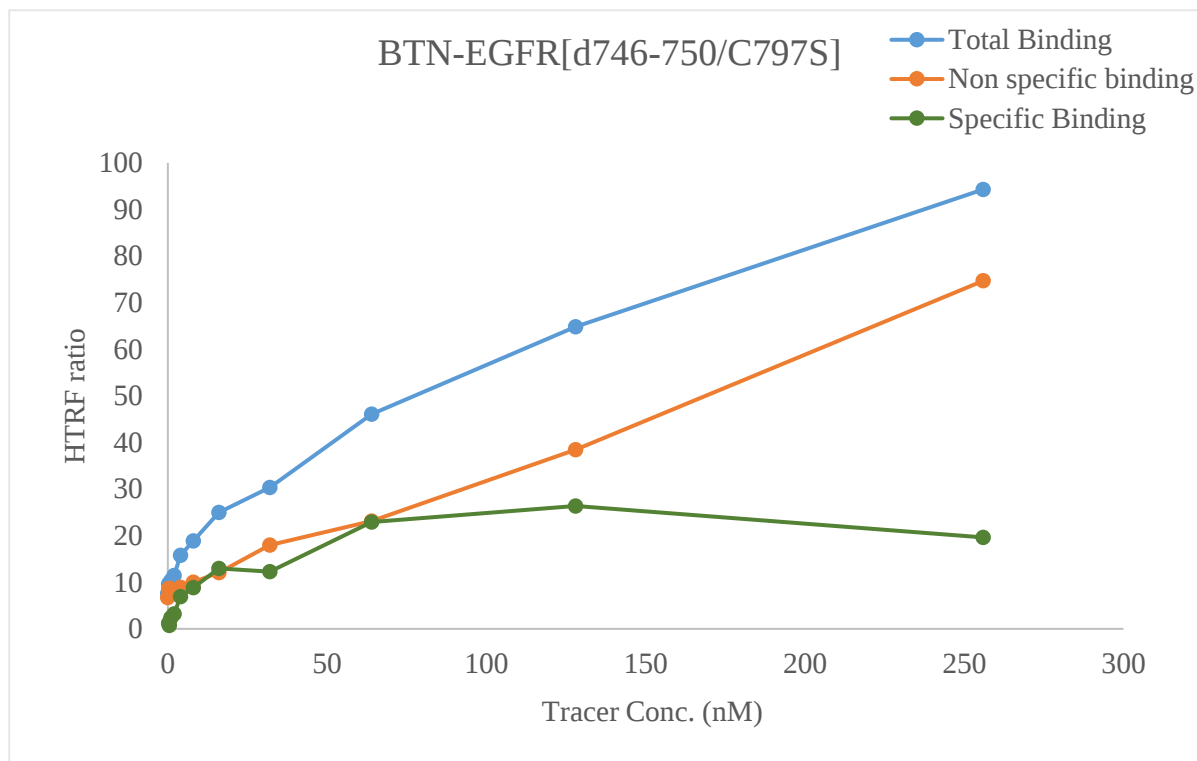


Kinase **BTN-EGFR[d746-750/C797S]**
 Cat.# 08-764-20N
 Tracer revvity Inc.
 Dasatinib-Red (#62KB02REDC)

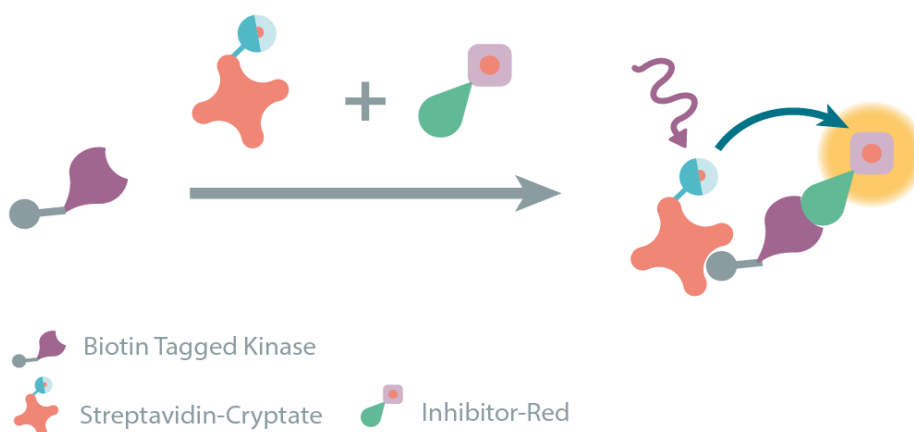


KD (nM) 14

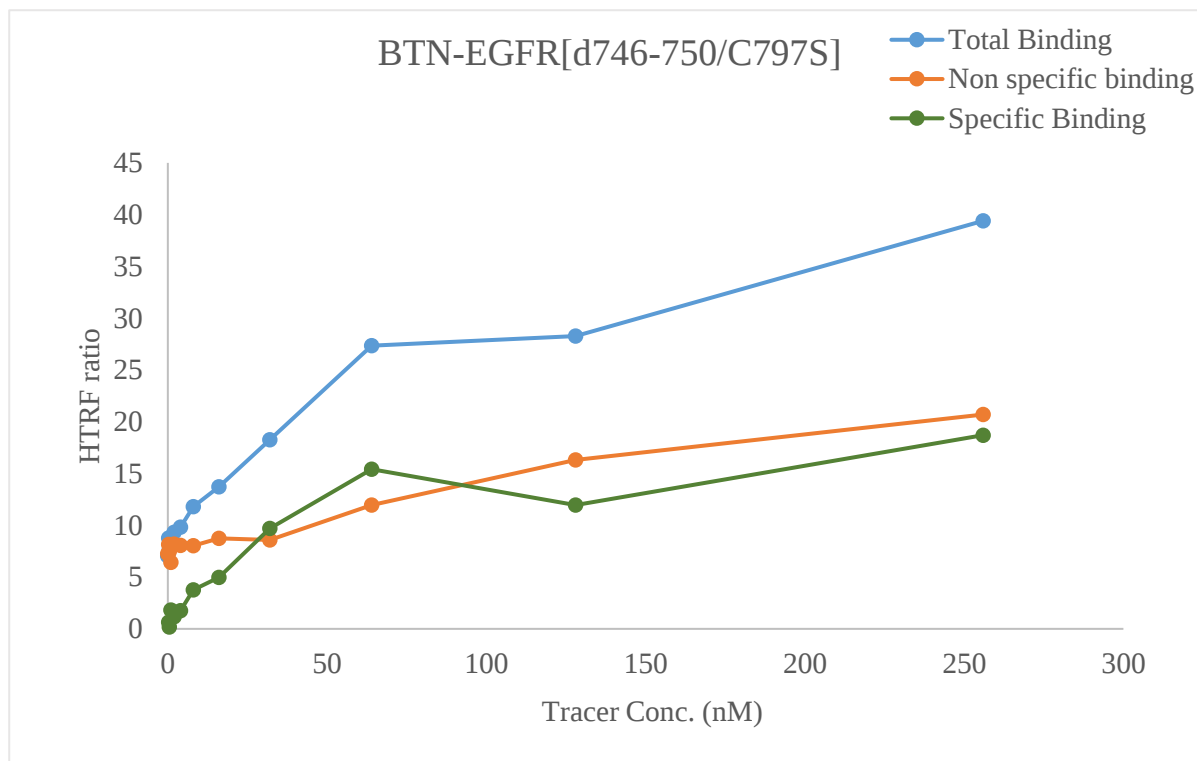
Kinase conc. 5 nM
 Eu labeled-streptavidin 5 nM

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[d746-750/C797S]**
 Cat.# 08-764-20N
 Tracer revvity Inc.
 Staurosporine-Red (#62KB01REDC)



KD (nM) 35

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

