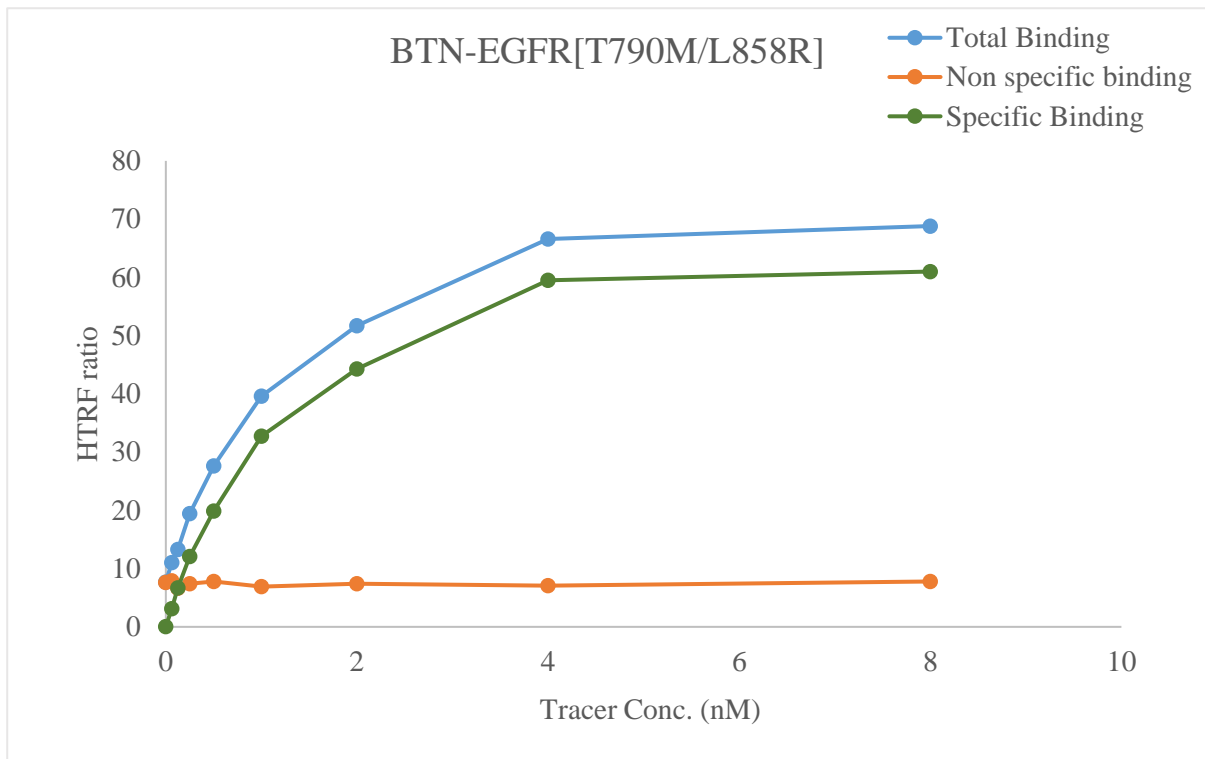


Kinase **BTN-EGFR[T790M/L858R]**

Cat.# 08-410-20N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



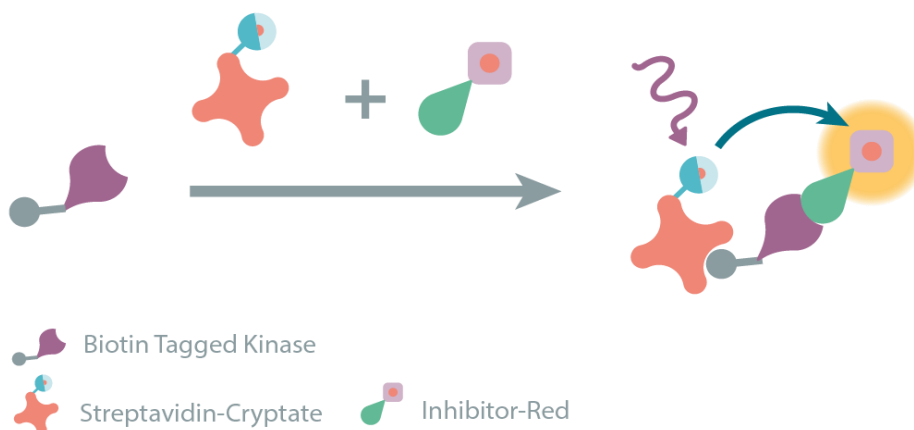
KD (nM) 1.1

Kinase conc. 5 nM

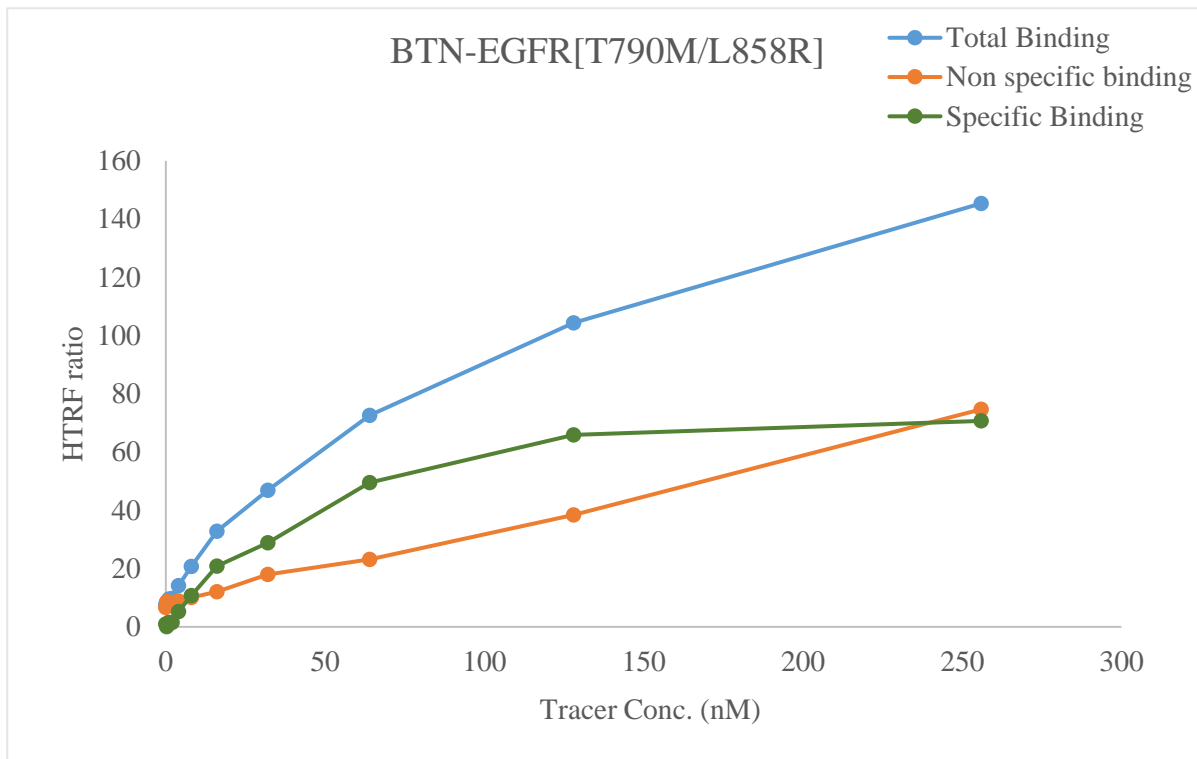
Eu labeled-streptavidin 5 nM

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[T790M/L858R]**
 Cat.# 08-410-20N
 Tracer revvity Inc.
 Dasatinib-Red (#62KB02REDC)



KD (nM) 55

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

