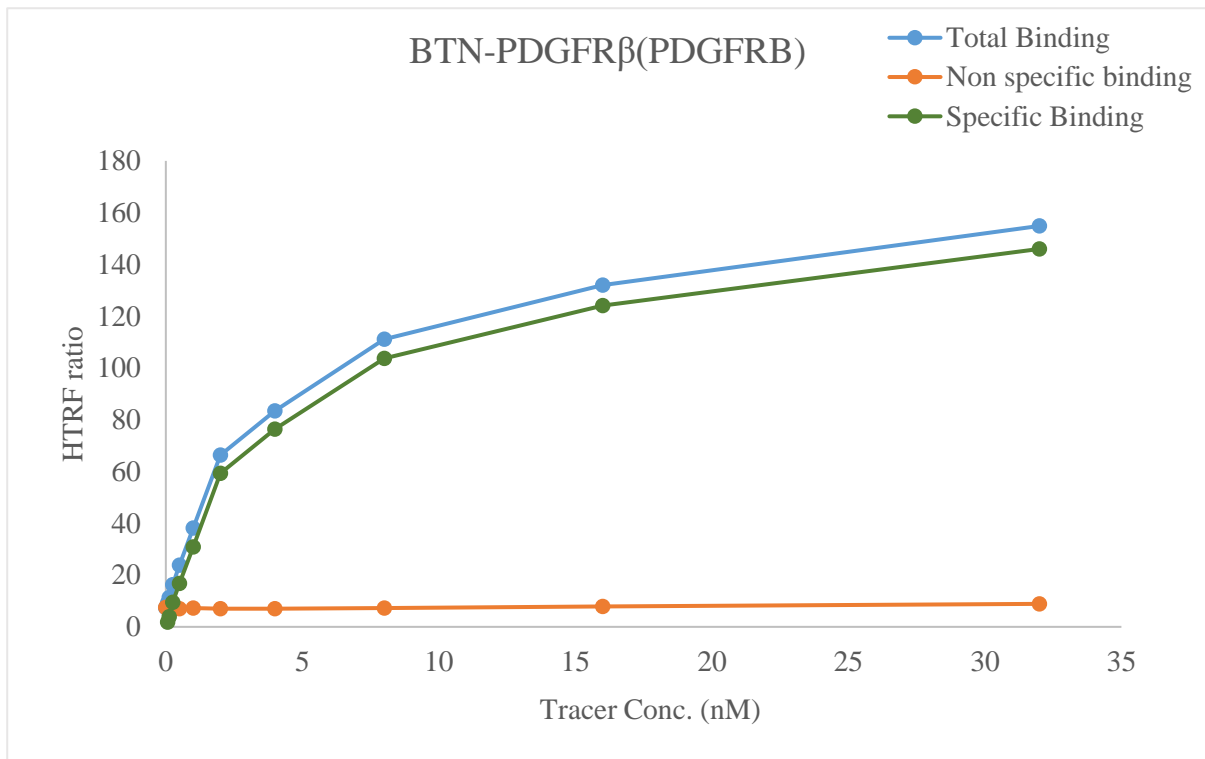


Kinase **BTN-PDGFR β (PDGFRB)**

Cat.# 08-458-20N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



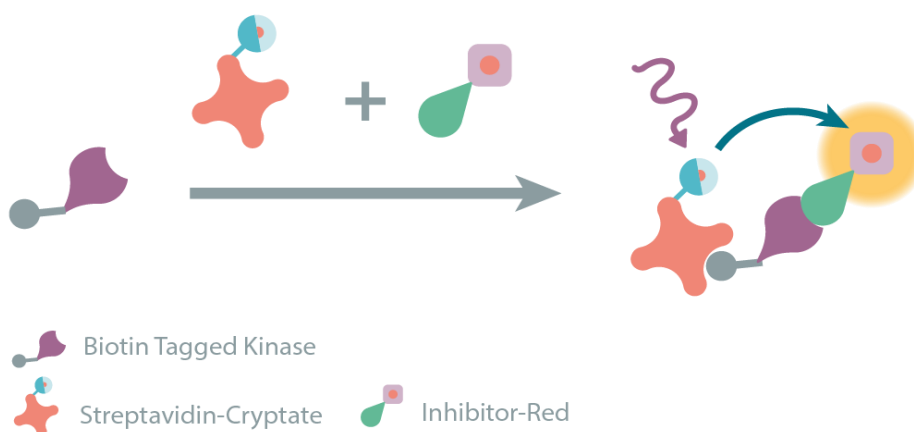
KD (nM) 4.6

Kinase conc. 5 nM

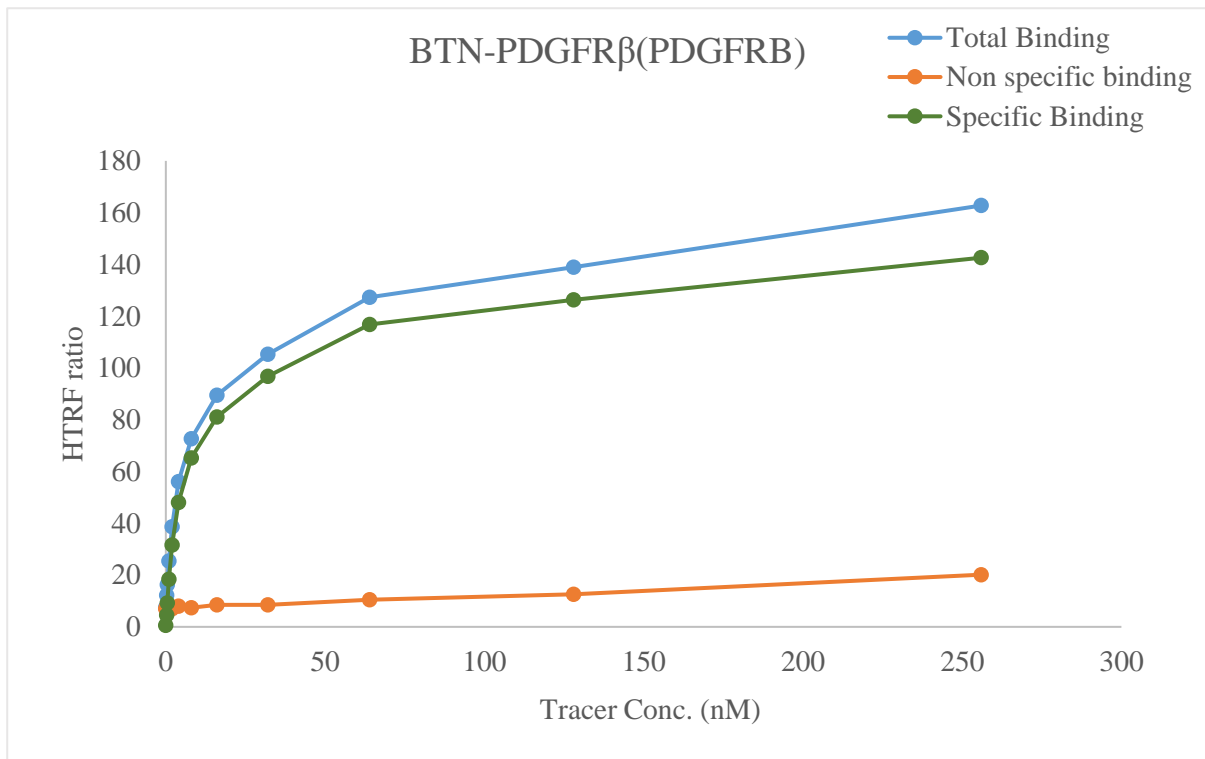
Eu labeled-streptavidin 5 nM

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-PDGFR β (PDGFRB)**
 Cat.# 08-458-20N
 Tracer revvity Inc.
 Sunitinib-Red (#62KB03REDC)



KD (nM) 8.2

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

