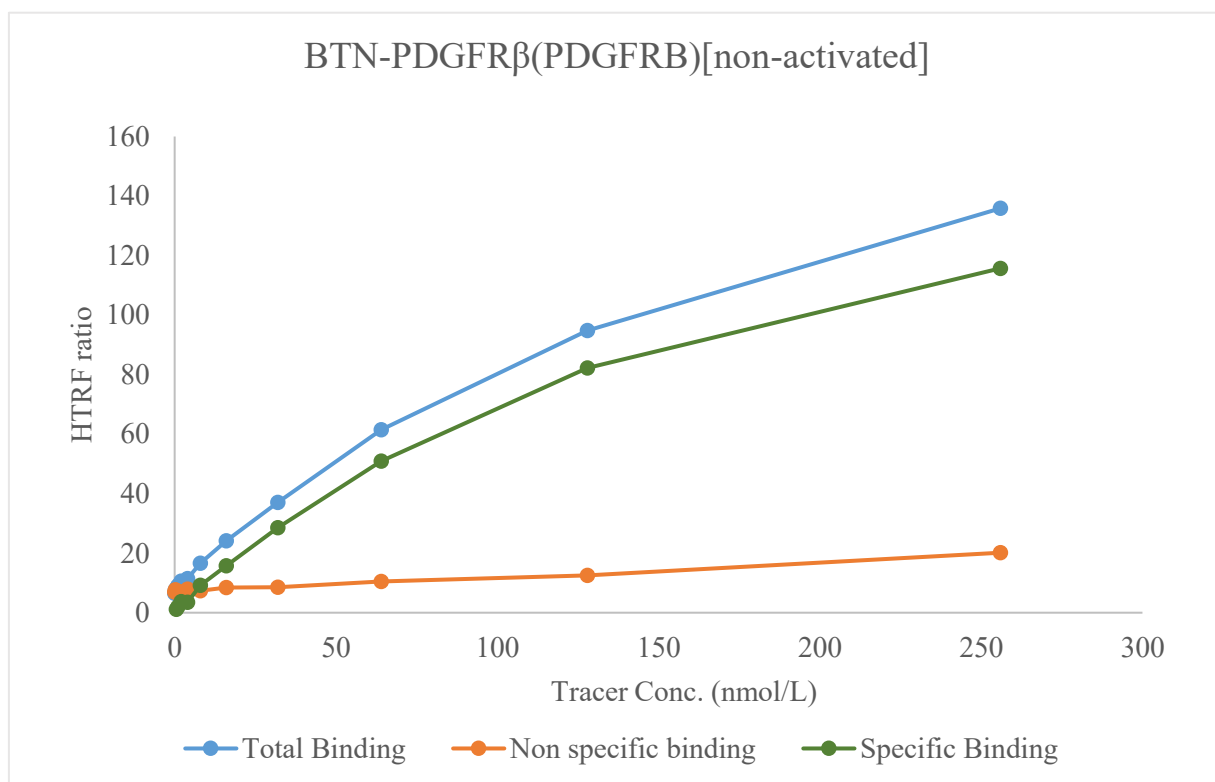


Kinase **BTN-PDGFR β (PDGFRB)[non-activated]**

Cat.# 08-458-23N



Tracer	revvity Inc. Sunitinib-Red, (#62KB03REDC)
KD	190 nmol/L
Kinase Conc.	5 nmol/L
Eu labeled-streptavidin	5 nmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	-
Reaction Time	1 Hour