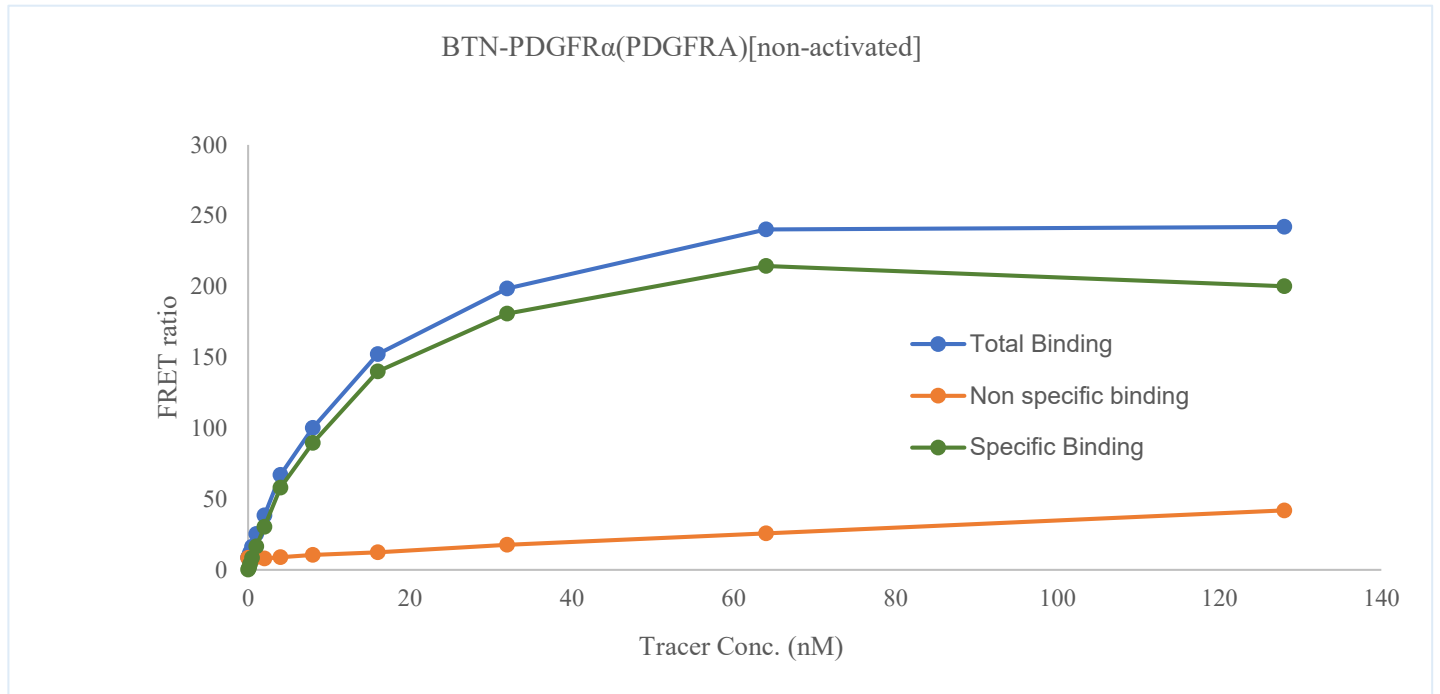


Kinase **BTN-PDGFR α (PDGFRA)[non-activated]**
 Cat.# 08-457-23N
 Tracer Revvity Inc.
 Dasatinib-Red (#62KB02REDC)



KD (nM)	12.1
Kinase conc.	5nM
Eu labeled-streptavidin	5nM

Follow the guide to set up your HTRF® binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF® kinase binding

