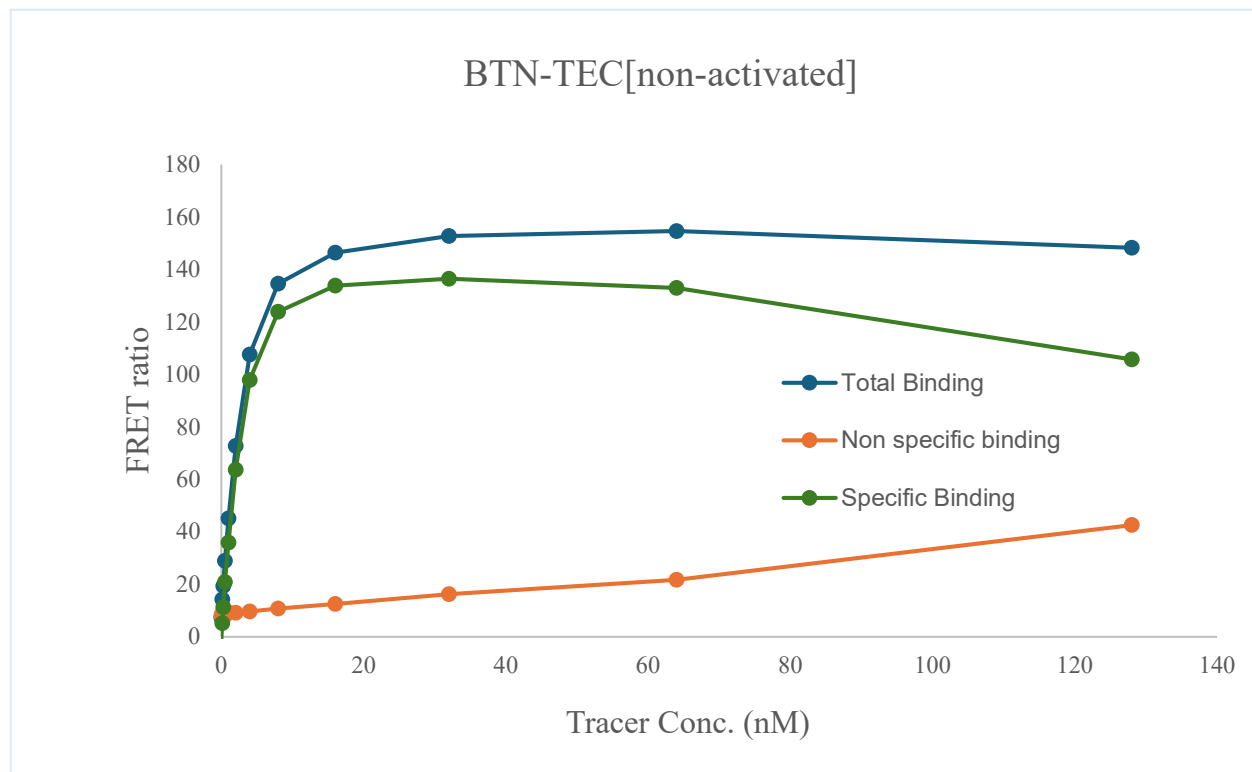


Kinase **BTN-TEC[non-activated]**
 Cat.# 08-482-23N
 Tracer Revvity Inc.
 Dasatinib-Red (#62KB02REDC)



KD (nM) 2.22
 Kinase conc. 5 nM
 Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF® binding assay

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

