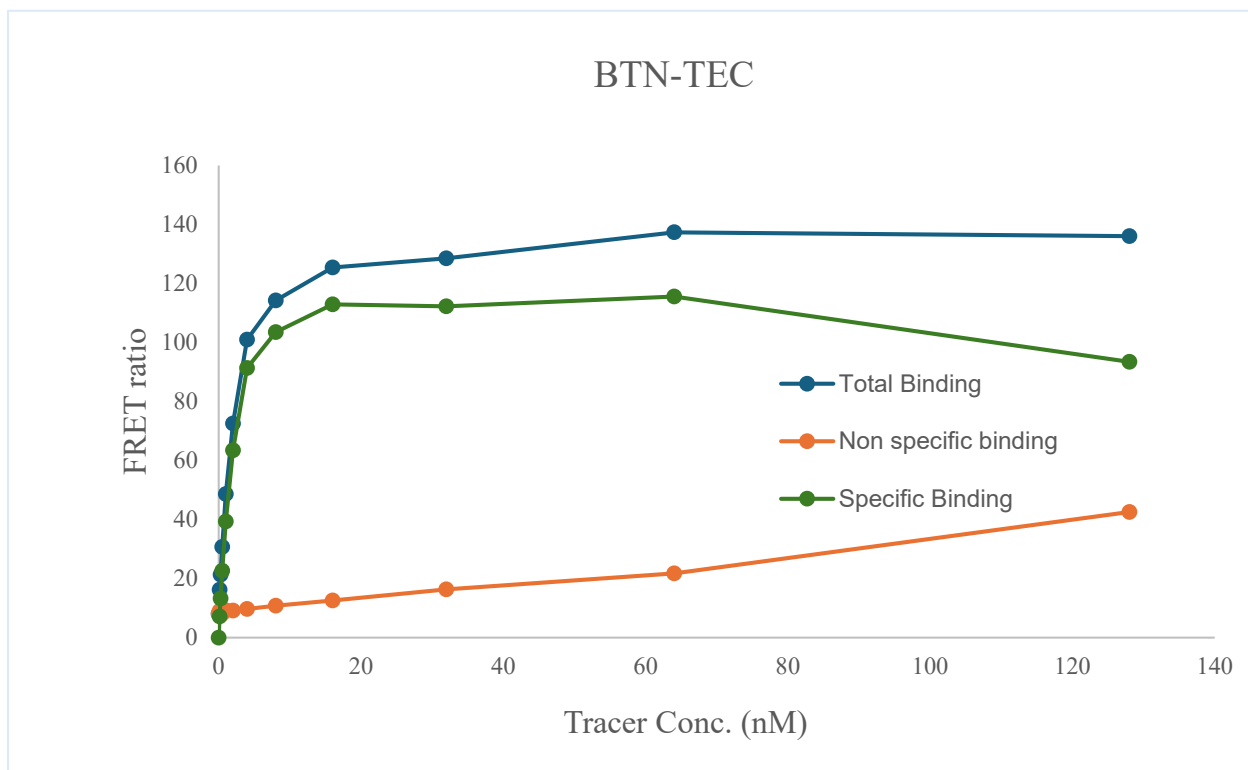


Kinase **BTN-TEC**
 Cat.# 08-482-20N
 Tracer Revvity Inc.
 Dasatinib-Red (#62KB02REDC)



KD (nM) 1.69
 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

