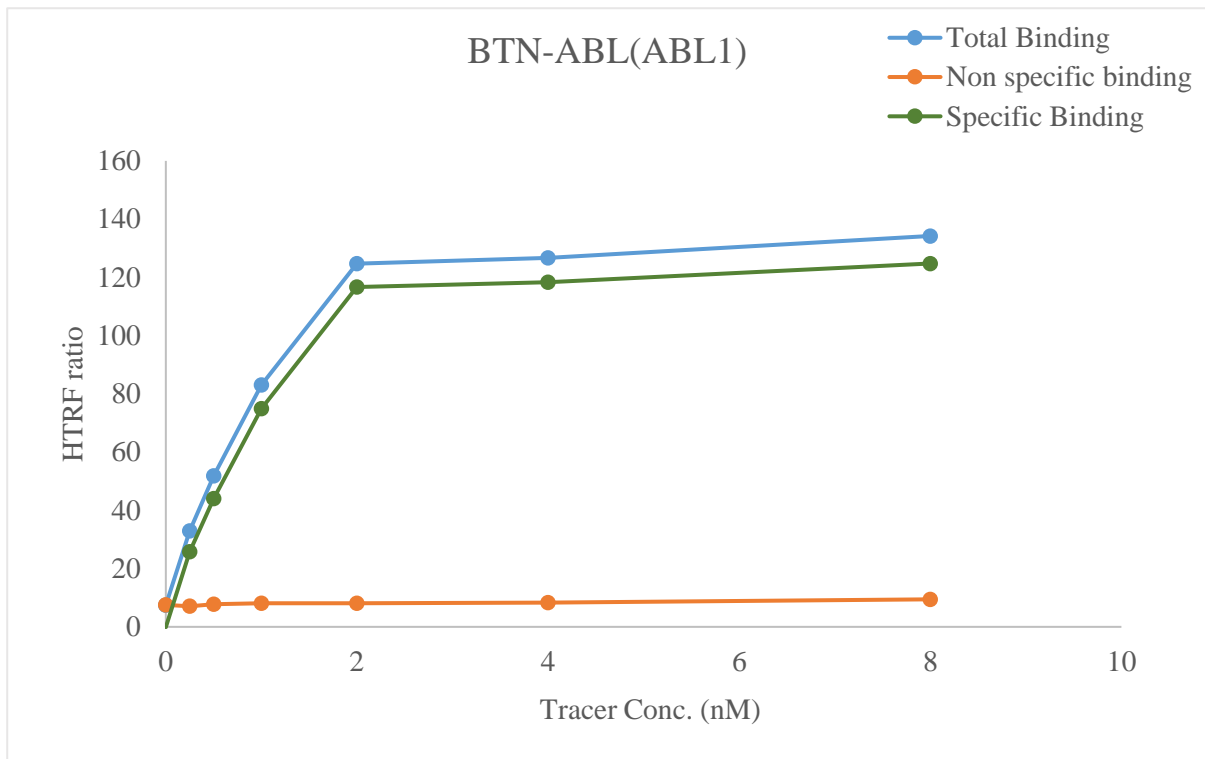


Kinase **BTN-ABL(ABL1)**
 Cat.# 08-401-20N
 Tracer revvity Inc.
 Dasatinib-Red (#62KB02REDC)



KD (nM) 0.76

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

