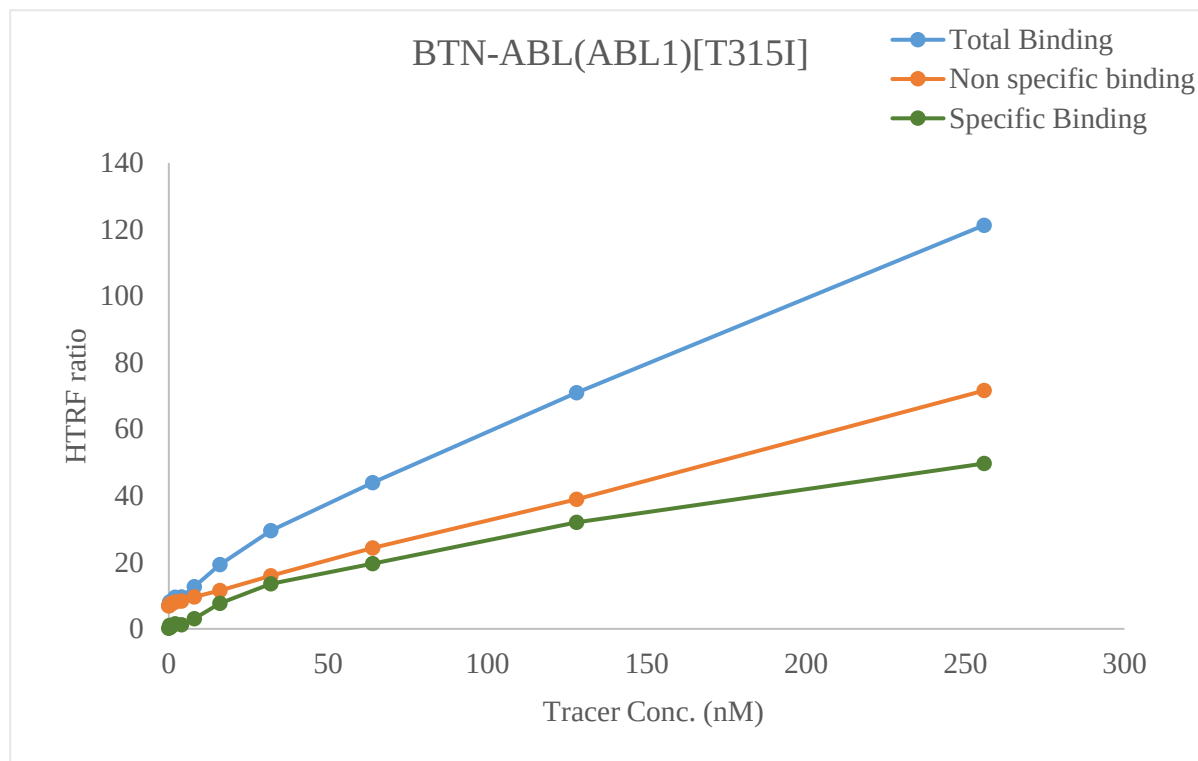


Kinase **BTN-ABL(ABL1)[T315I]**

Cat.# 08-493-20N

Tracer revvity Inc.

Dasatinib-Red (#62KB02REDC)



KD (nM) 210

Kinase conc. 5 nM

Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF<sup>®</sup> binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF<sup>®</sup> kinase binding

