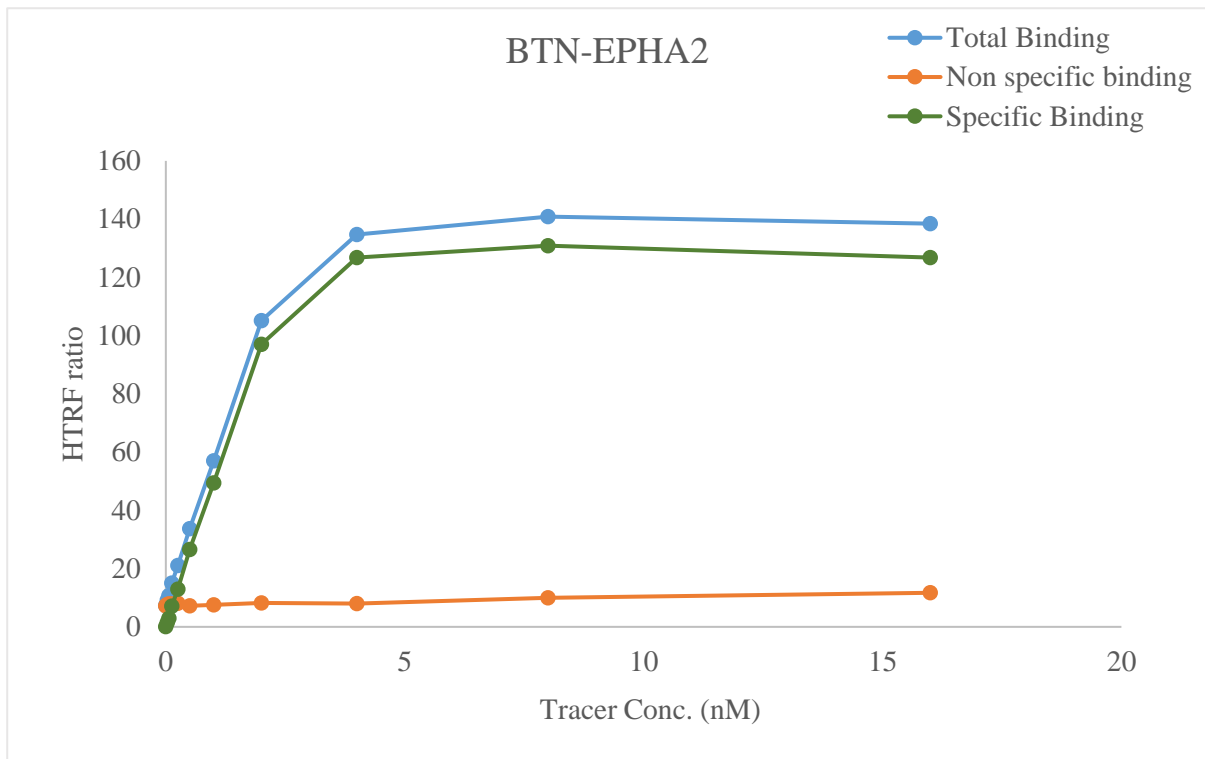


Kinase **BTN-EPHA2**

Cat.# 08-421-20N

Tracer revvity Inc.

Dasatinib-Red (#62KB02REDC)



KD (nM) 1.5

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

