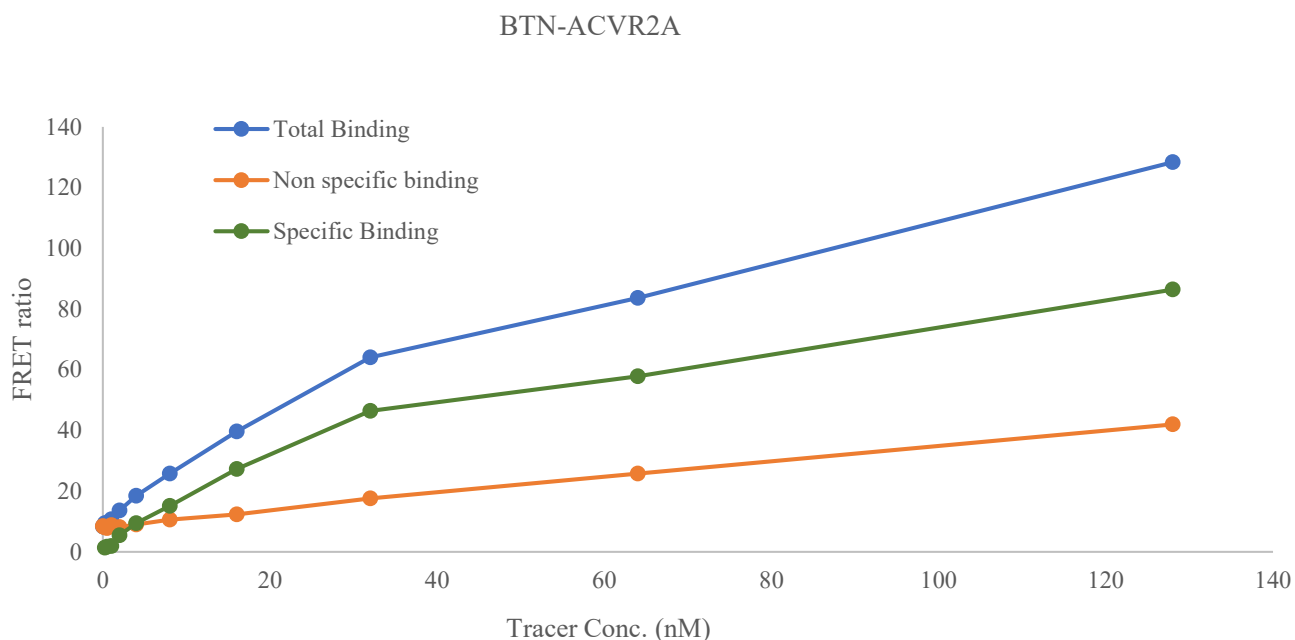


Kinase **BTN-ACVR2A**
 Cat.# 09-431-20N
 Tracer Revvity Inc.
 Dasatinib-Red (#62KB02REDC)



KD (nM)	55.2
Kinase conc.	5nM
Eu labeled-streptavidin	5nM

Follow the guide to set up your HTRF® binding assay

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

