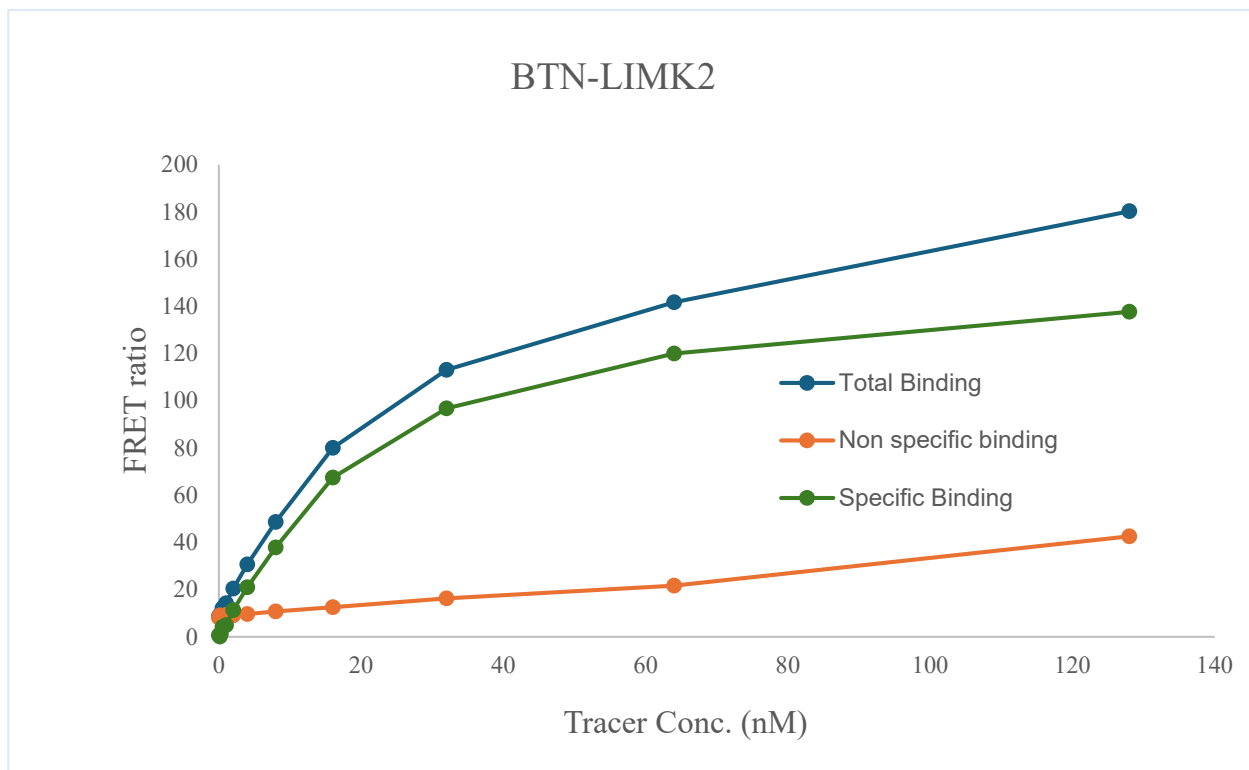


Kinase **BTN-LIMK2**  
 Cat.# 09-406-20N  
 Tracer Revvity Inc.  
 Dasatinib-Red (#62KB02REDC)



KD (nM) 24.3  
 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

