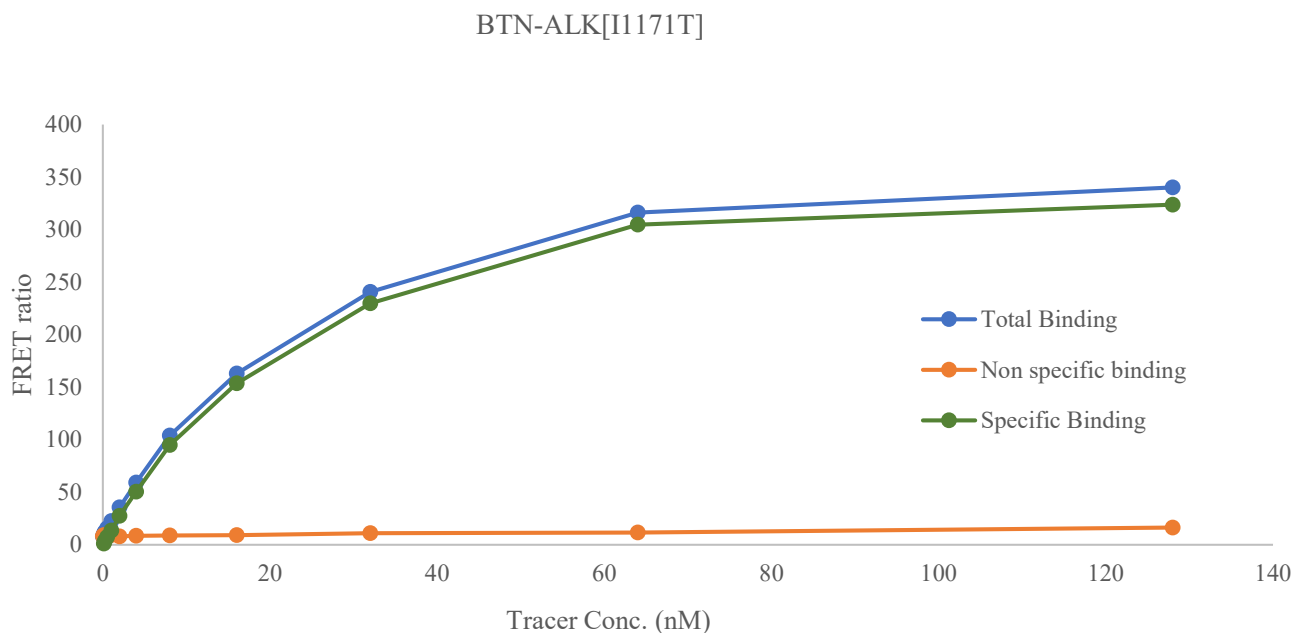


Kinase **BTN-ALK[I1171T]**
 Cat.# 08-790-20N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB01REDC)



KD (nM)	24.9
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

