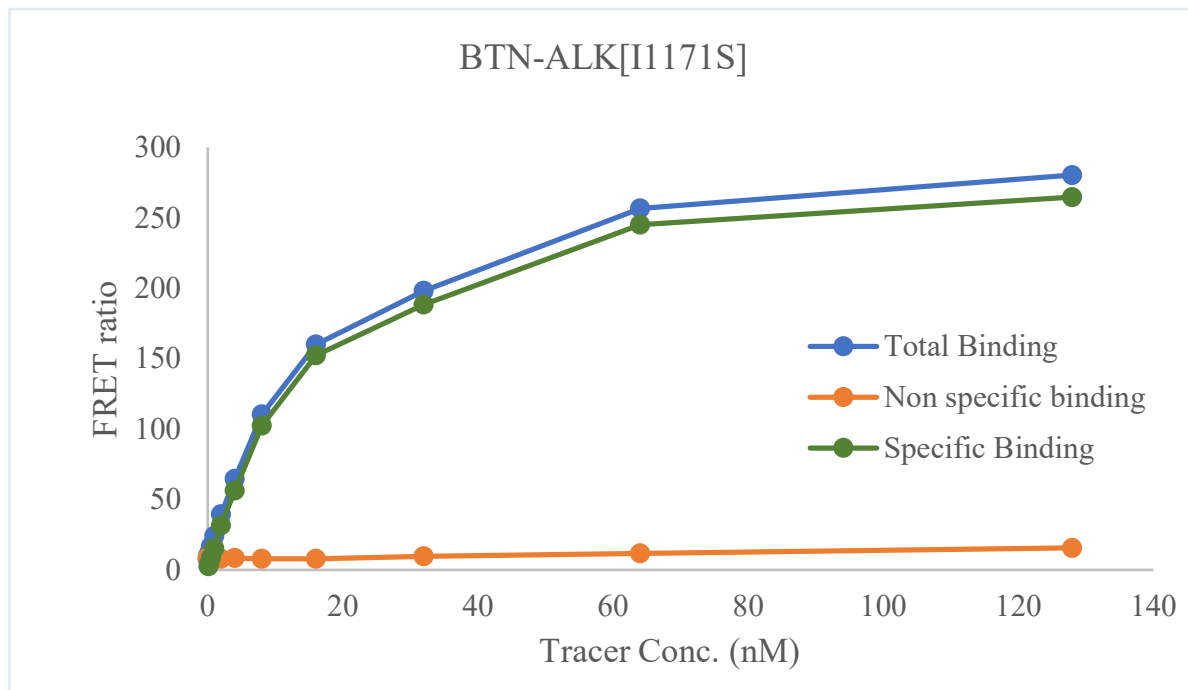


Kinase **BTN-ALK[I1171S]**

Cat.# 08-769-20N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM) 16.51

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

