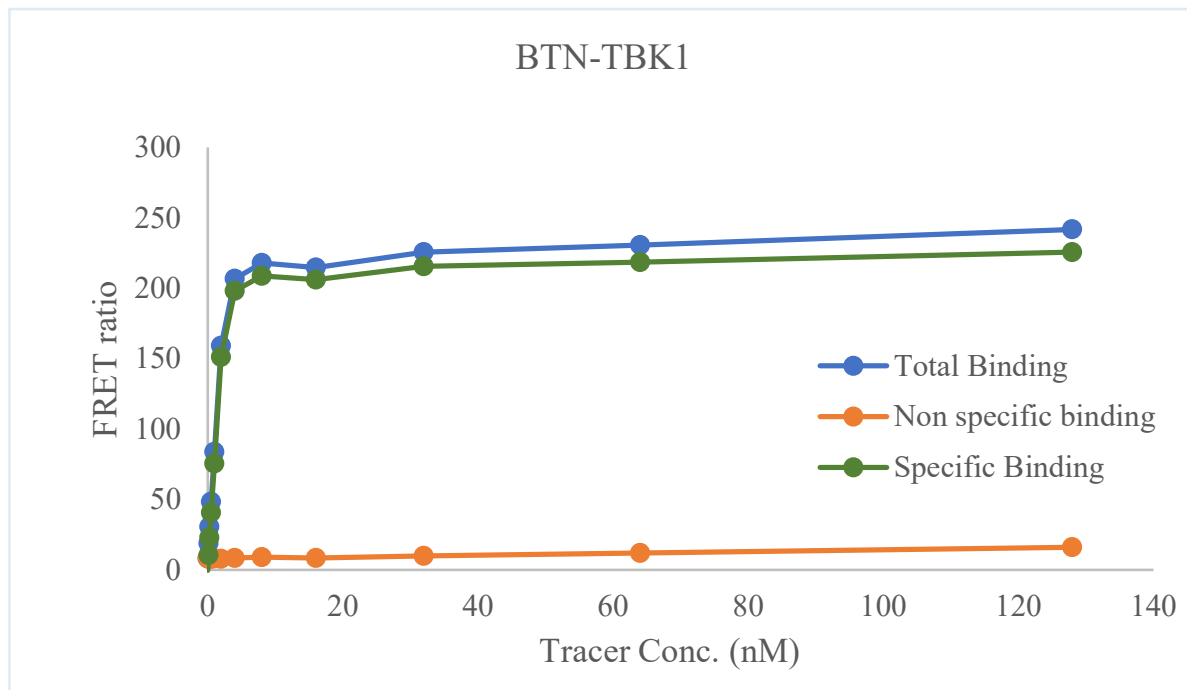


Kinase **BTN-TBK1**
 Cat.# 05-415-20N
 Tracer Revvity Inc.
 Staurosporine-Red



KD (nM) 1.47
 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

