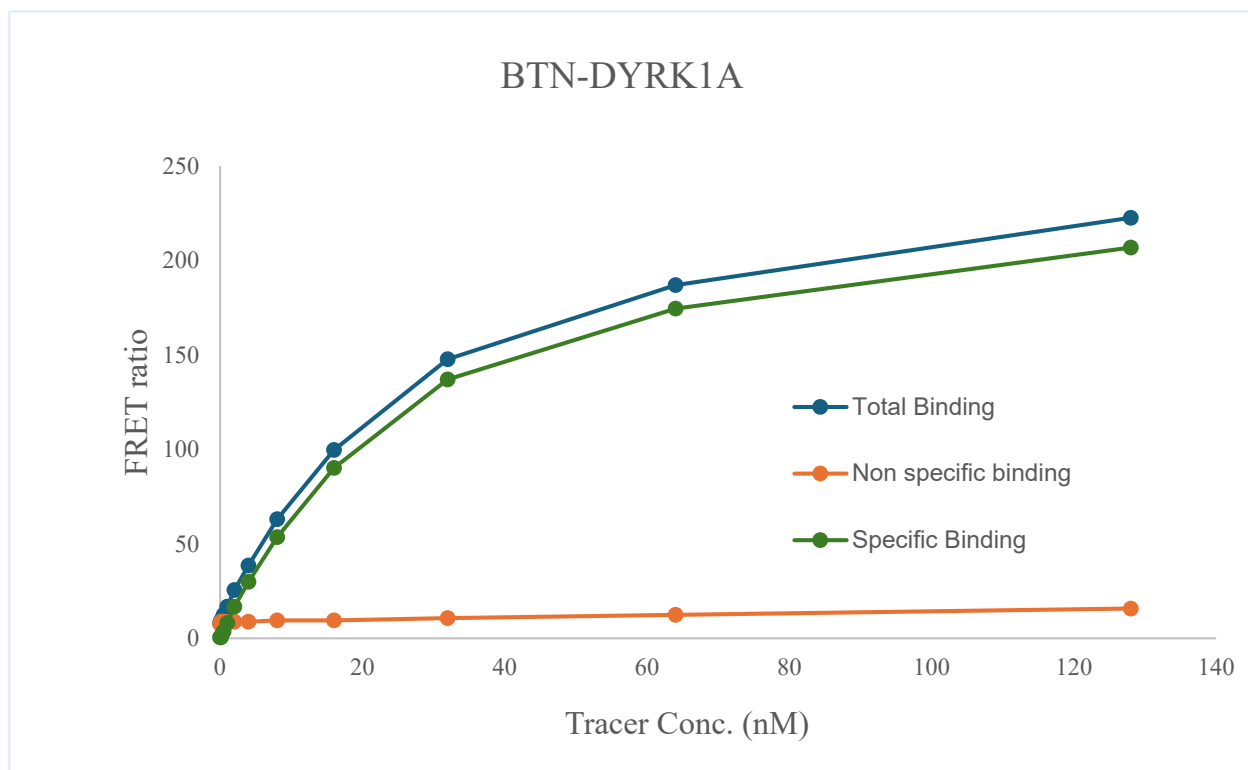


Kinase **BTN-DYRK1A**
 Cat.# 04-430-20N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB02REDC)



KD (nM) 28.4
 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

