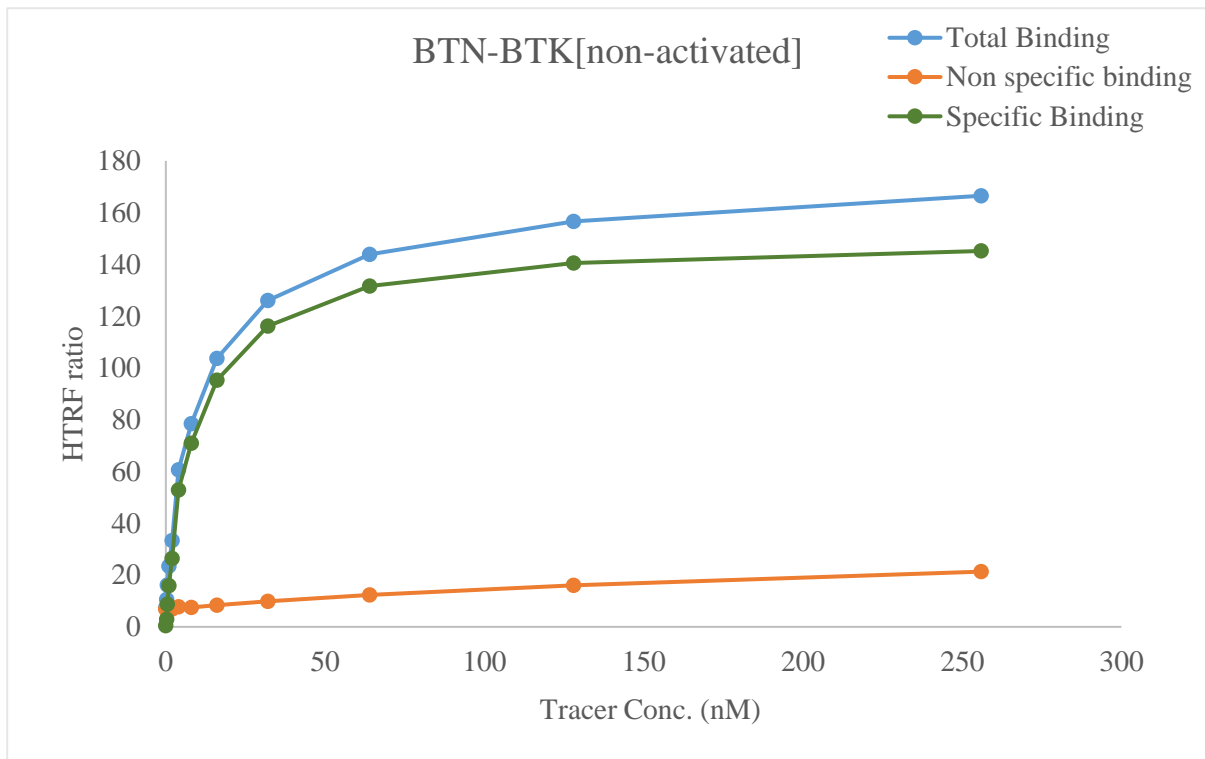


Kinase **BTN-BTK[non-activated]**

Cat.# 08-480-23N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 8.6

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

