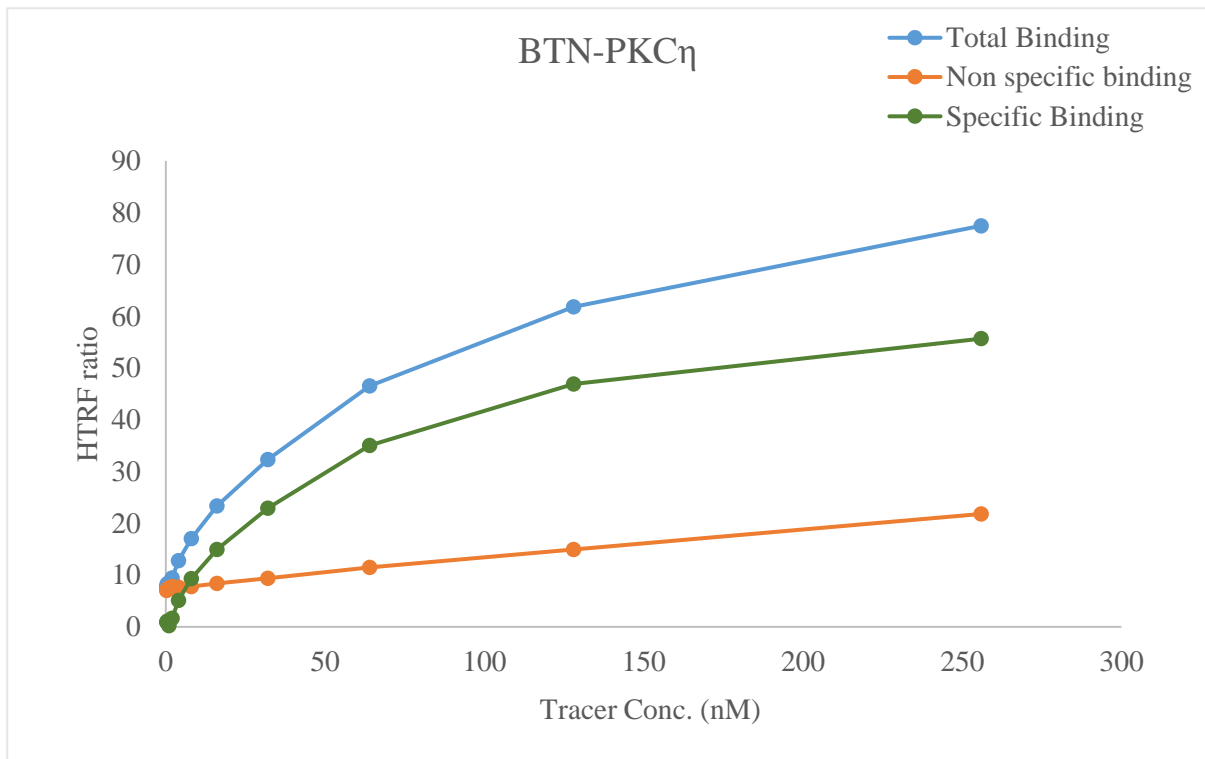


Kinase **BTN-PKC $\eta$**

Cat.# 01-438-20N

Tracer revvity Inc.

Staurosporine-Red (#62KB01REDC)



KD (nM) 59

Kinase conc. 5 nM

Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF<sup>®</sup> binding assay

Determination of the dissociation constant (Kd) of the fluorescent tracer using HTRF<sup>®</sup> kinase binding

