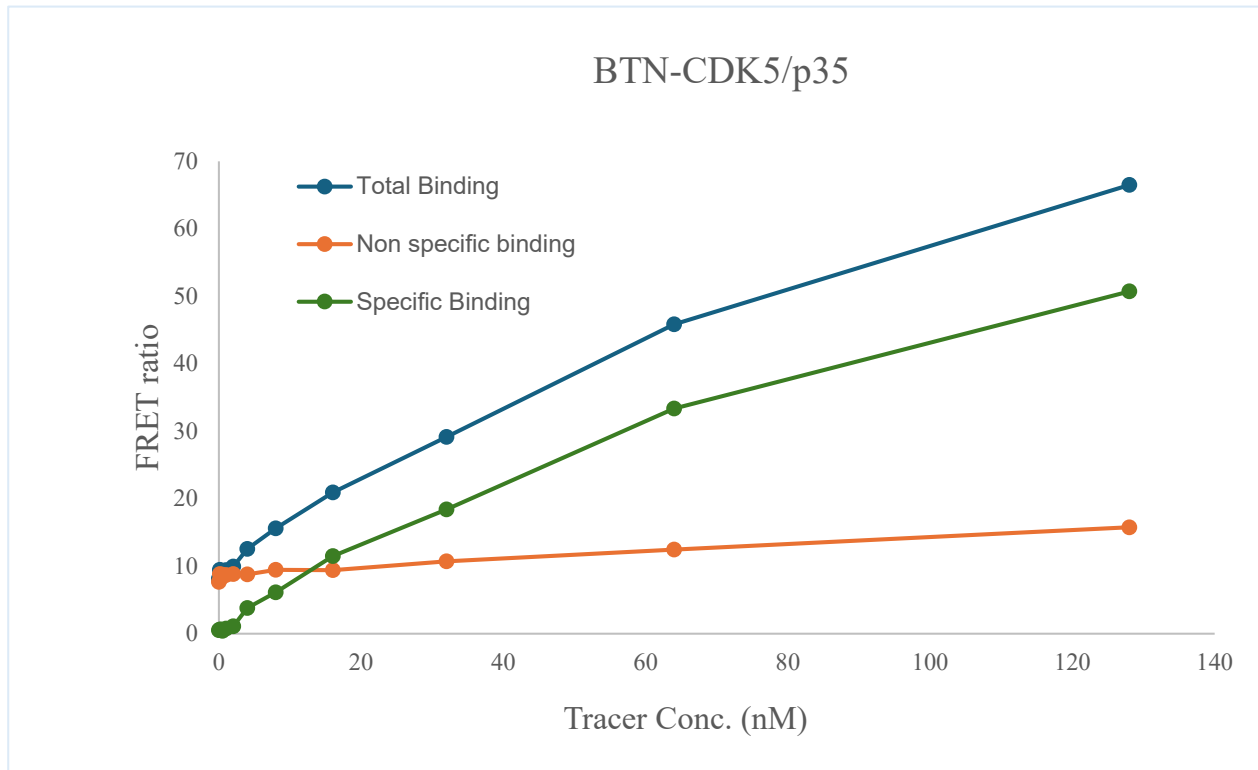


Kinase **BTN-CDK5/p35**
 Cat.# 04-492-20N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB02REDC)



KD (nM)	140
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

