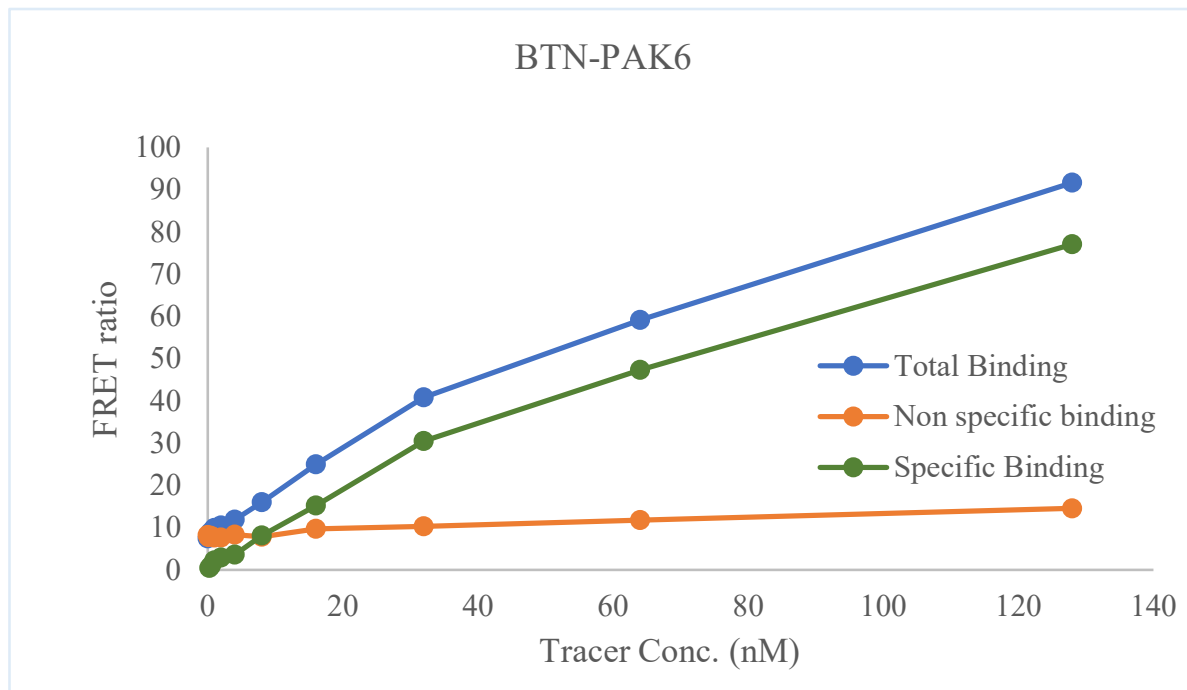


Kinase **BTN-PAK6**
 Cat.# 07-428-20N
 Tracer Revvity Inc.
 Staurosporine-Red



KD (nM)	151
Kinase conc.	5nM
Eu labeled-streptavidin	5nM

Follow the guide to set up your HTRF® binding assay

Determination of the dissociation constant (K_d) of the fluorescent tracer using HTRF® kinase binding

