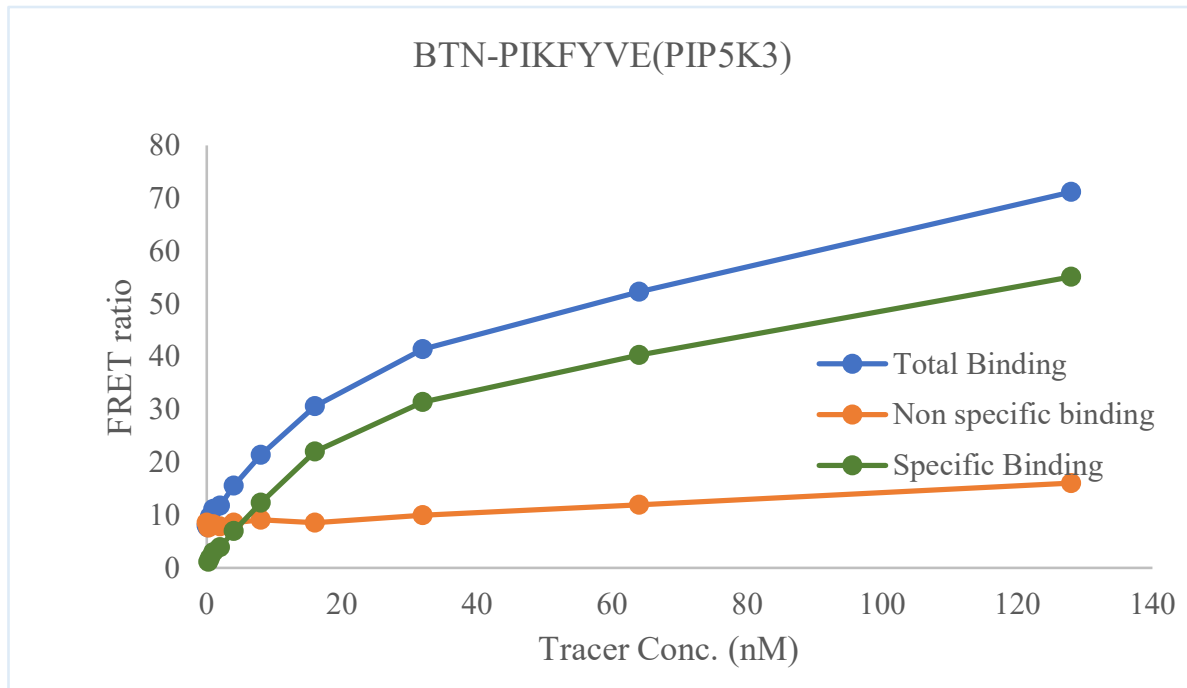


Kinase **BTN-PIKFYVE(PIP5K3)**

Cat.# 11-418-20N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM)	38.3
Kinase conc.	5nM
Eu labeled-streptavidin	5nM

Follow the guide to set up your HTRF® binding assay



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

