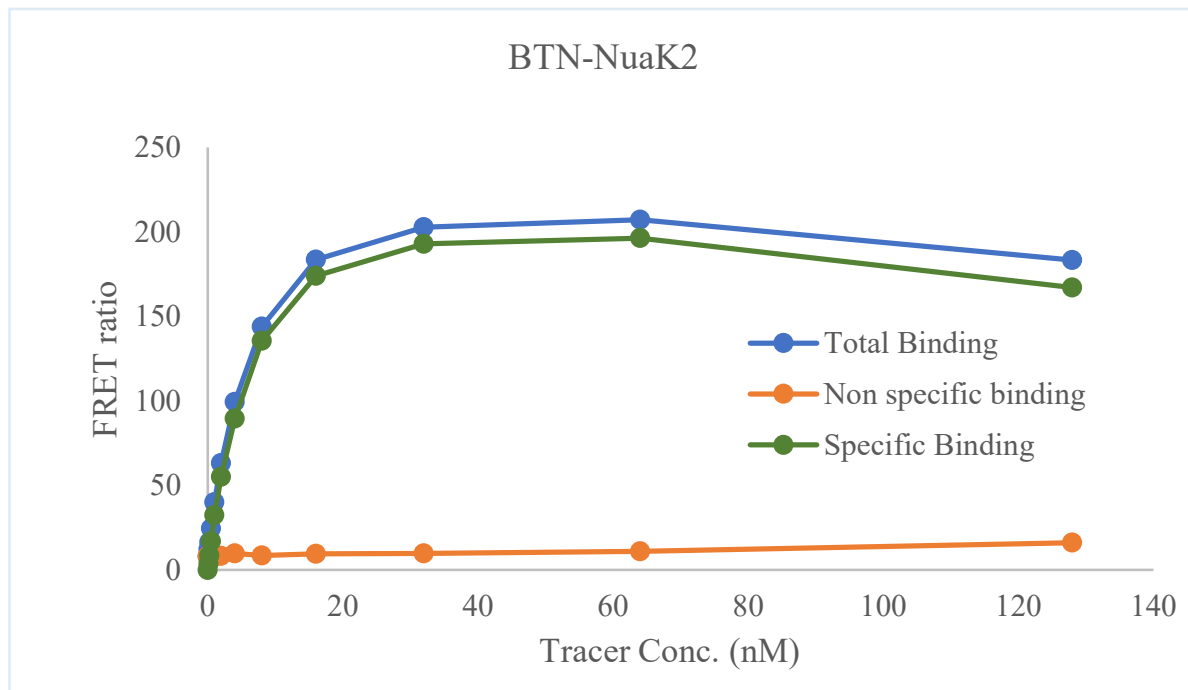


Kinase **BTN-NuaK2**
 Cat.# 02-427-20N
 Tracer Revvity Inc.
 Staurosporine-Red



KD (nM)	4.78
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

