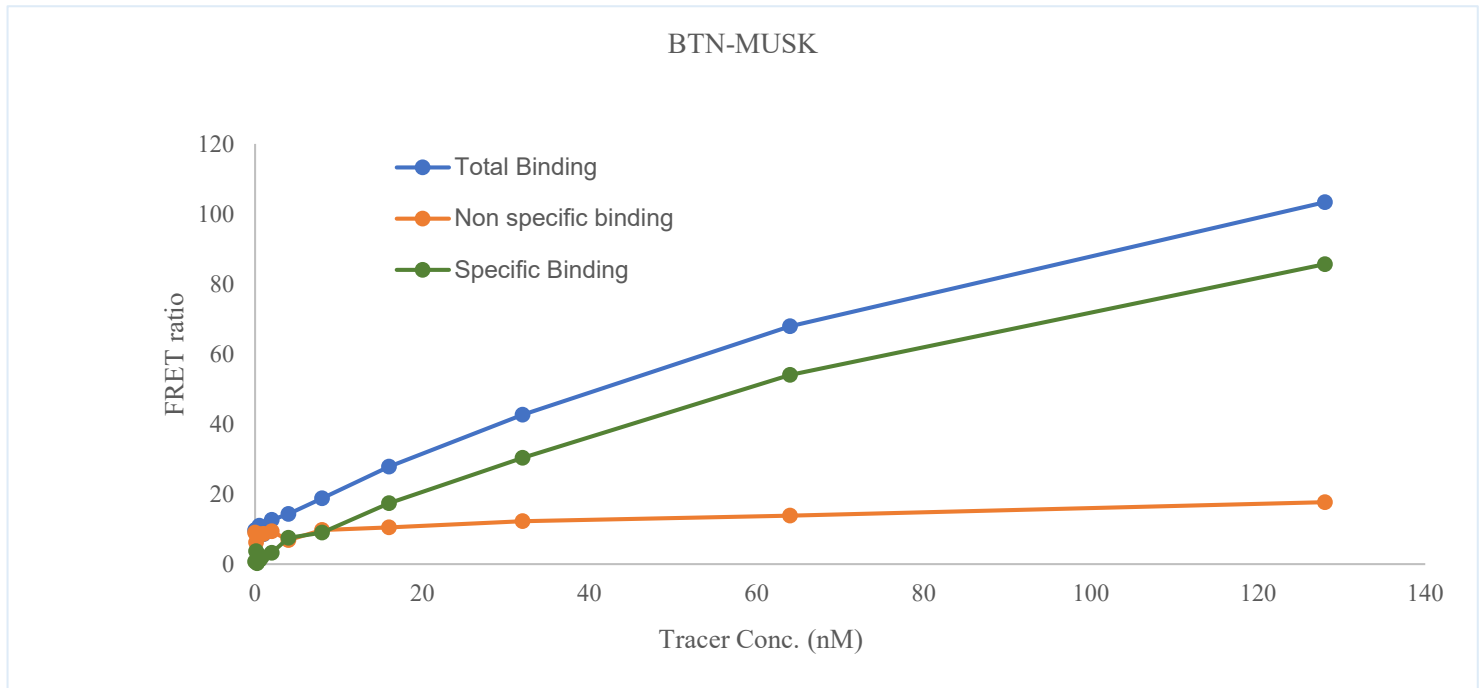


Kinase **BTN-MUSK**
 Cat.# 08-453-20N
 Tracer Revvity Inc.
 Staurosporine-Red (#62KB01REDC)



	KD (nM)	159
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

