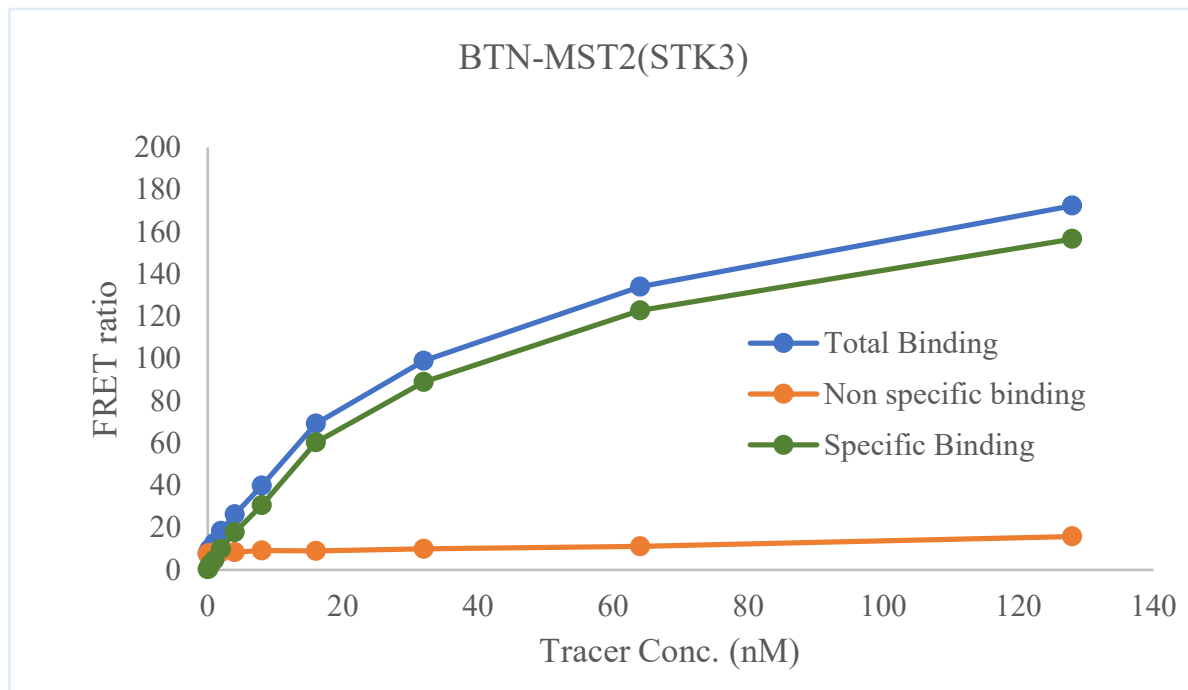


Kinase **BTN-MST2(STK3)**

Cat.# 07-417-20N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM) 41.5

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

