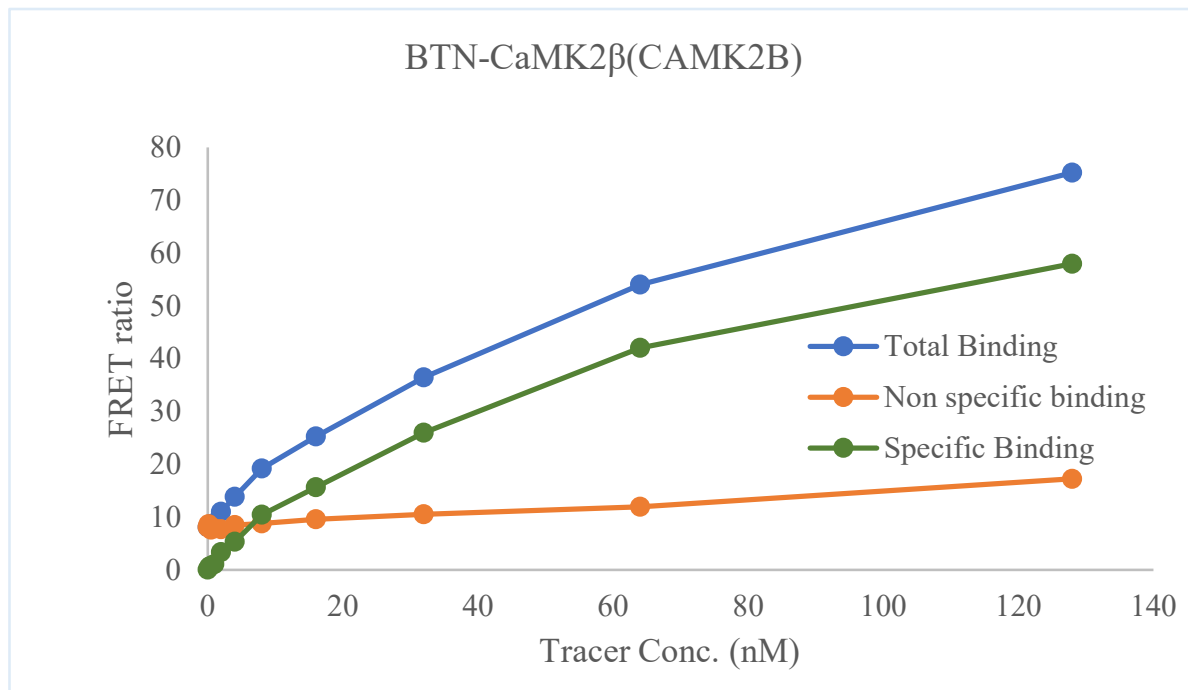


Kinase **BTN-CaMK2 β (CAMK2B)**

Cat.# 02-410-20N

Tracer Revvity Inc.

Staurosporine-Red



KD (nM) 77.9

Kinase conc. 5 nM

Eu labeled-streptavidin 5 nM

Follow the guide to set up your HTRF® binding assay



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

