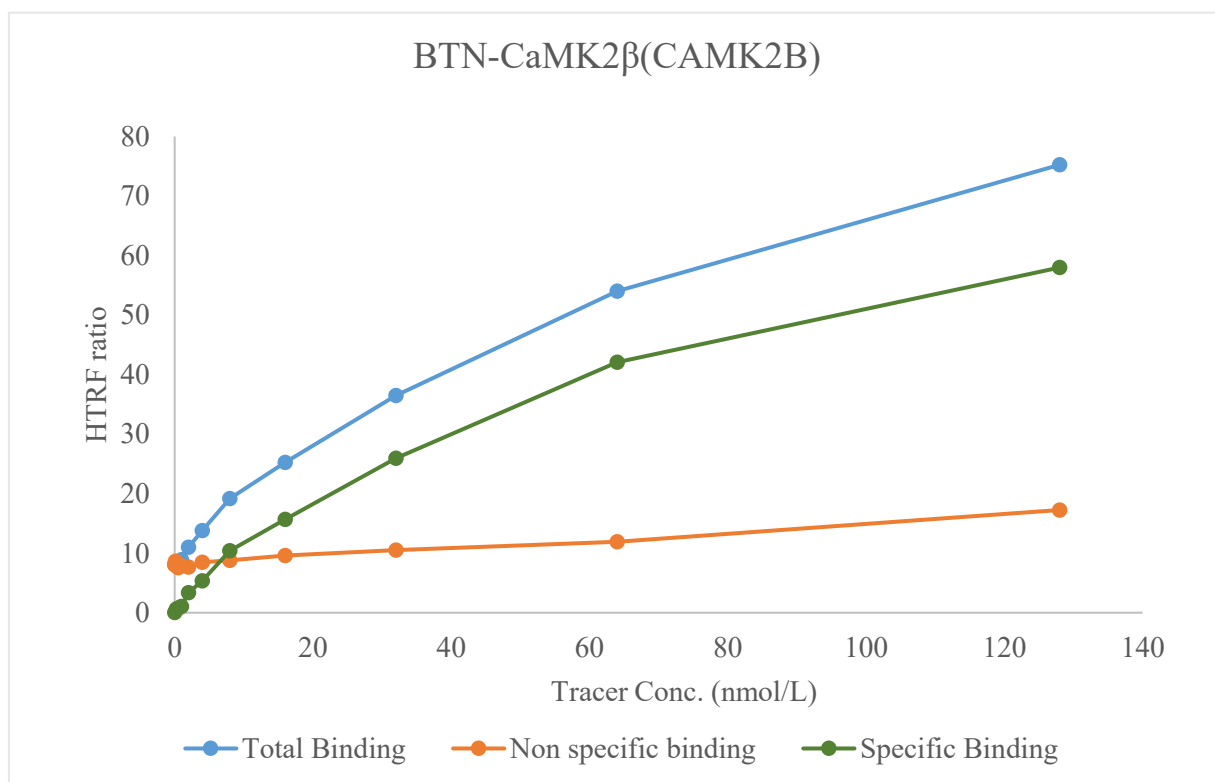


Kinase **BTN-CaMK2 β (CAMK2B)**

Cat.# 02-410-20N



Tracer	revvity Inc. Staurosporine-Red, (#62KB01REDC)
KD	78 nmol/L
Kinase Conc.	5 nmol/L
Eu labeled-streptavidin	5 nmol/L
Metal	Mg
Metal Conc.	5 mmol/L
Additives	Calmodulin 0.01 mg/mL
Reaction Time	1 Hour