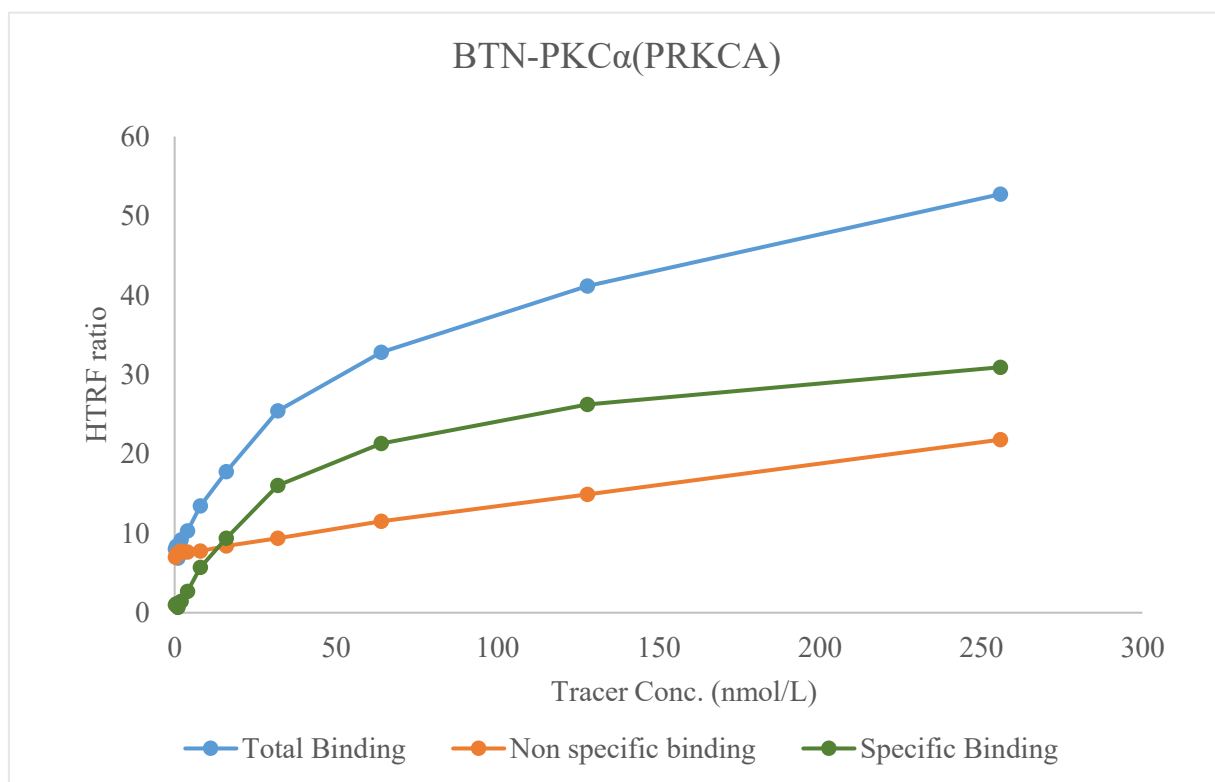


Kinase **BTN-PKC α (PRKCA)**

Cat.# 01-433-20N



Tracer	revvity Inc. Staurosporine-Red, (#62KB01REDC)
KD	41 nmol/L
Kinase Conc.	5 nmol/L
Eu labeled-streptavidin	5 nmol/L
Metal	Mg, Ca
Metal Conc.	5, 0.05 mmol/L
Additives	5 µg/mL Diacyl Glycerol, 50 µg/mL Phosphatid
Reaction Time	1 Hour