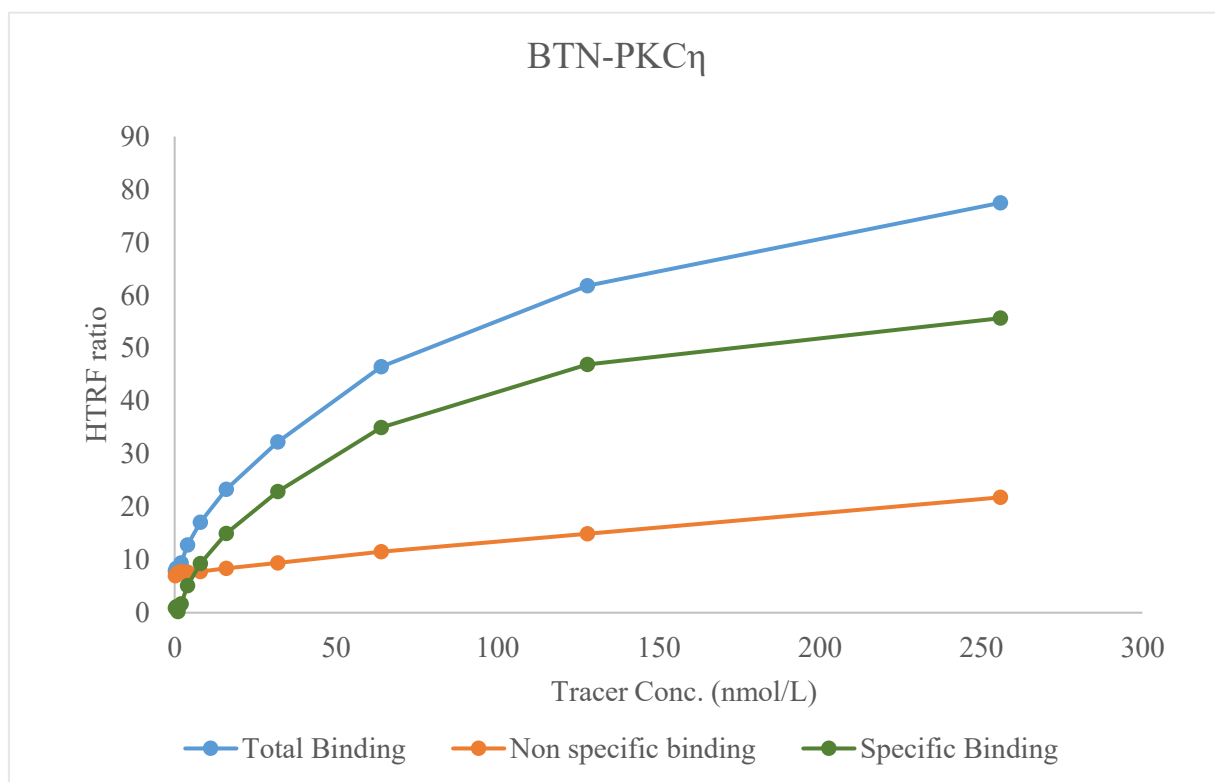


Kinase **BTN-PKC $\eta$**

Cat.# 01-438-20N



|                         |                                                  |
|-------------------------|--------------------------------------------------|
| Tracer                  | revvity Inc.<br>Staurosporine-Red, (#62KB01REDC) |
| KD                      | 59 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                                           |