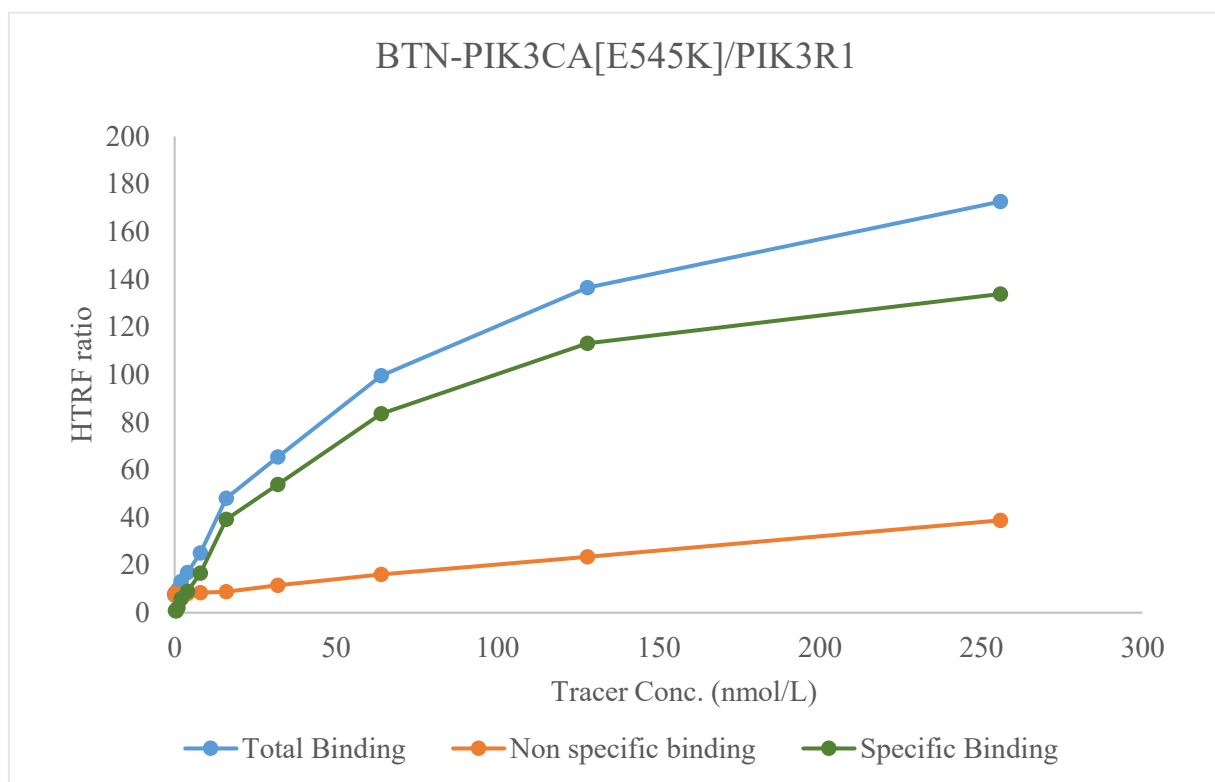


Kinase **BTN-PIK3CA[E545K]/PIK3R1**

Cat.# 11-414-20N



Tracer

Thermo Scientific

Kinase Tracer 1710, PV6088

KD

62 nmol/L

Kinase Conc.

5 nmol/L

Eu labeled-streptavidin

5 nmol/L

Metal

Mg

Metal Conc.

5 mmol/L

Additives

-

Reaction Time

1 Hour