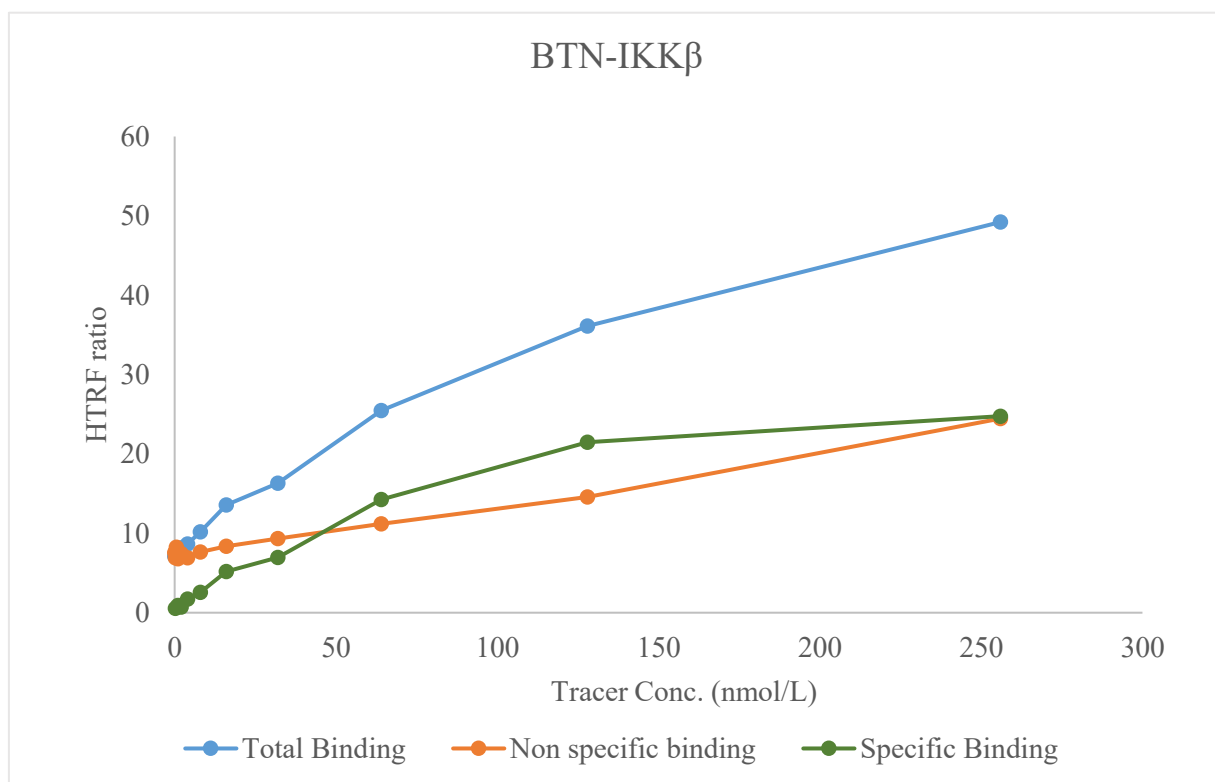


Kinase **BTN-IKK β**

Cat.# 05-484-20N



Tracer

Thermo Scientific

Kinase Tracer 236, PV5592

KD

110 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