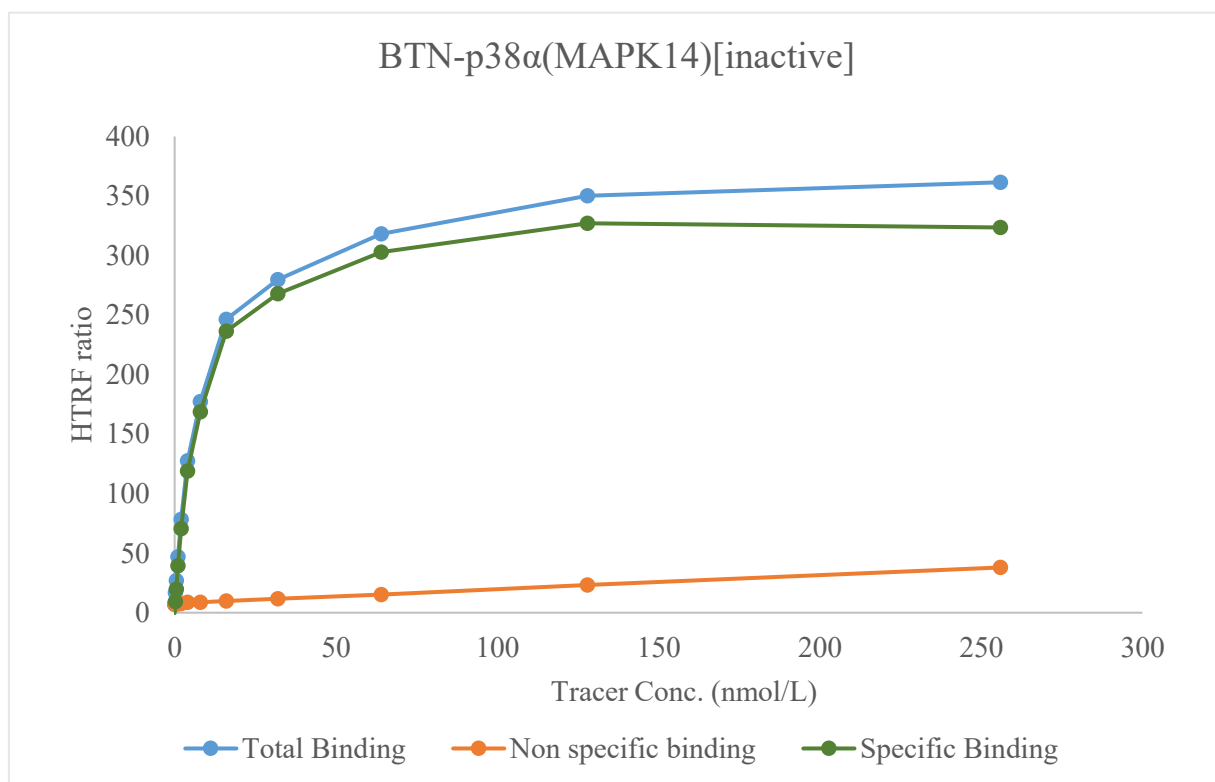


Kinase **BTN-p38 α (MAPK14)[inactive]**

Cat.# 04-452-10-20N



Tracer

Thermo Scientific

Kinase Tracer 199, PV5830

KD

7.6 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