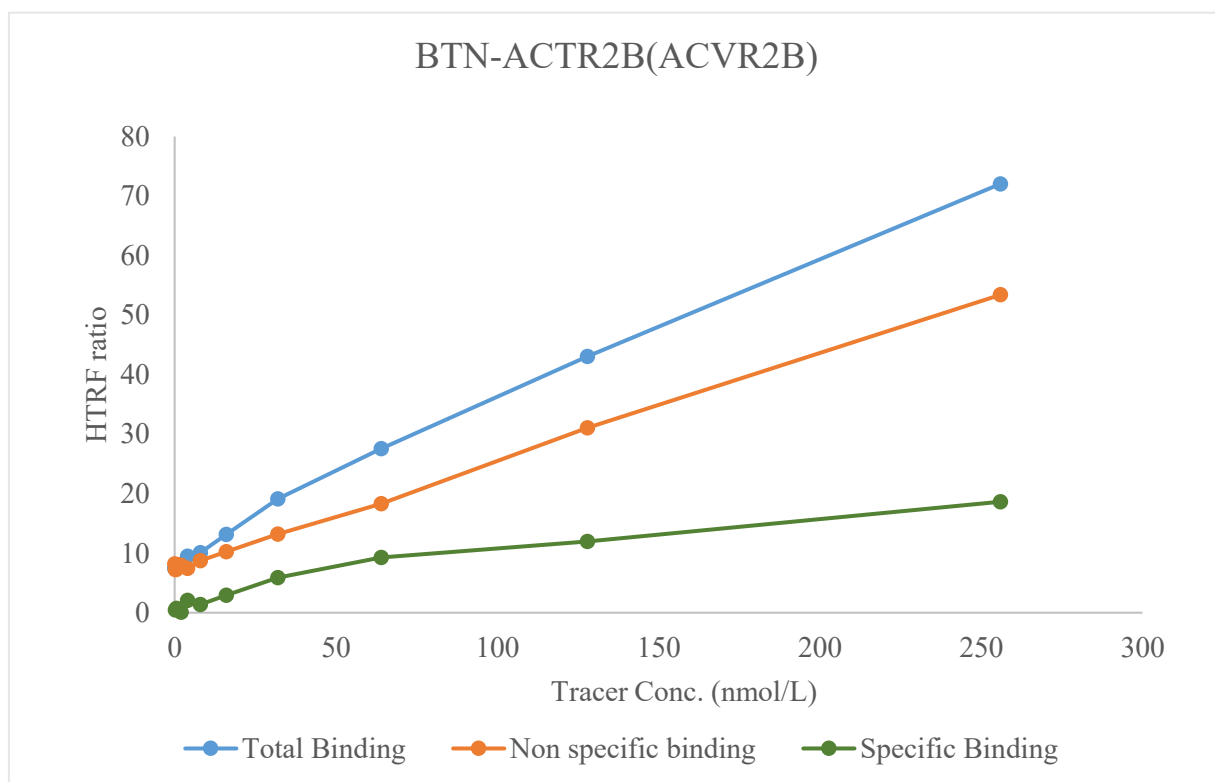


Kinase **BTN-ACTR2B(ACVR2B)**

Cat.# 09-432-20N



Tracer

Thermo Scientific

Kinase Tracer 178, PV5593

KD

150 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